

PRODUCT DATASHEET

ChemiScreen™ GABA_B GABA Family Receptor Stable Cell Line

CATALOG NUMBER: HTS119C

CONTENTS: 2 vials of mycoplasma-free cells, 1 ml per vial.

STORAGE: Vials are to be stored in liquid N₂.

BACKGROUND

ChemiScreen cell lines are constructed in the Chem-1 host, which supports high levels of functional receptor expression on the cell surface. Chem-1 cells contain high endogenous levels of Gα15, a promiscuous G protein, allowing most receptors to couple to the calcium signaling pathway.

The neurotransmitter γ -aminobutyric acid (GABA) exerts its effects through an ion channel, GABA_A, and a GPCR, GABA_B. Functional GABA_B is a heterodimer composed of the GABA_{B1} and GABA_{B2} subunits, which share 35% sequence identity and belong to the class 3 family of GPCRs. The GABA_{B1} subunit, which exists as splice variants GABA_{B1a} and GABA_{B1b}, binds directly to GABA and is required for agonist activation. The GABA_{B2} and GABA_{B1} subunits associate by formation of a coiled coil by their C-terminal tails; this association masks an ER retention sequence in GABA_{B1} to permit export from the ER and trafficking to the cell surface. In addition to its chaperone function, GABA_{B2} is the component that couples to G_i to reduce intracellular cAMP. Agonists of GABA_B, such as baclofen, are used clinically for treatment of muscle spasticity, migraine headache and musculoskeletal pain (Bowery *et al.*, 2002). Cloned human GABA_{B1b}/GABA_{B2} receptor-expressing ChemiScreen cells were constructed by stable transfection of Chem-1 cells with GABA_{B1b}/GABA_{B2} which supports high levels of recombinant GABA_{B1b}/GABA_{B2} expression on the cell surface and contains high levels of the promiscuous G protein Gα15 to couple the receptor to the calcium signaling pathway. These stability-tested cells are ready for fluorescence-based assays for agonists, antagonists and modulators at the GABA_B receptor.

USE RESTRICTIONS

Please see **Limited Use Label License Agreement** (Label License Agreement) for further details.

WARNINGS

For Research Use Only; Not for Use in Diagnostic Procedures
Not for Animal or Human Consumption

GMO

This product contains genetically modified organisms.
Este producto contiene organismos genéticamente modificados.
Questo prodotto contiene degli organismi geneticamente modificati.
Dieses Produkt enthält genetisch modifizierte Organismen.
Ce produit contient organismes génétiquement des modifiés.
Dit product bevat genetisch gewijzigde organismen.
Tämä tuote sisältää geneettisesti muutettuja organismeja.
Denna produkt innehåller genetiskt ändrade organismer.

APPLICATIONS

Calcium Flux Fluorescence Assay

APPLICATION DATA

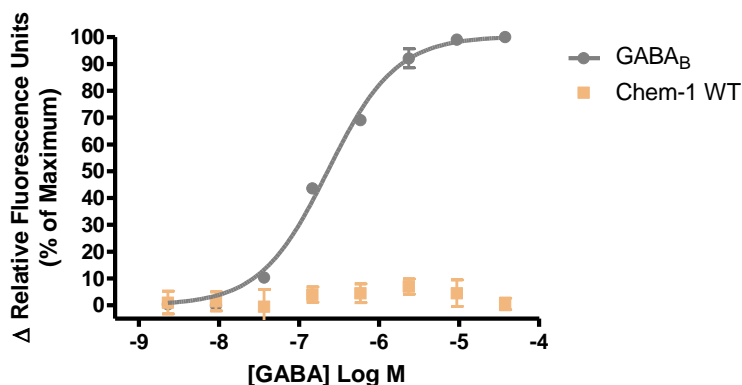


Figure 1. Representative data for activation of GABA_{B1b}/GABA_{B2} receptor stably expressed in Chem-1 cells induced by GABA using a fluorescent calcium flux assay. GABA_{B1b}/GABA_{B2}-expressing Chem-1 cells were seeded at 50,000 cells per well into a 96-well plate, and the following day the cells were loaded with a fluorescent calcium indicator. Calcium flux in response to the indicated ligand with a final concentration of 0.5% DMSO was determined on a Molecular Devices FLIPR^{TETRA}® with ICCD camera. Maximal fluorescence signal obtained in this experiment was 5,000 RLU. Similarly parental cells (catalog #: HTSCHEM-1) were tested to determine the specificity of the resulting signal.

Table 1. EC₅₀ values of GABA_{B1b}/GABA_{B2}-expressing Chem-1 cells.

LIGAND	ASSAY	POTENCY EC ₅₀ (nM)	REFERENCE
GABA	Calcium Flux - Fluorescence	230	Eurofins Internal Data

* The cell line was tested and found to have equivalent EC₅₀ and signal at 1, 3 and 6 weeks of continuous culture by calcium flux fluorescence. The Z' value, as defined with response to GABA, was 0.7.

CELL CULTURE

Table 2. Recommended Cell Culture Reagents (not provided)

Description	Component	Concentration	Supplier and Product Number
Basal Medium	DMEM high glucose Medium (4.5g/L)	-	Hyclone: SH30022
	Fetal Bovine Serum (FBS)	10%	Hyclone: SH3007003
	Non-Essential Amino Acids (NEAA)	1X	Hyclone: SH3023801
	HEPES	1X	EMD Millipore:TMS-003-C
Selection Medium	Basal Medium (see above)	-	
	Geneticin (G418)	250 µg/ml	Invivogen: ant-gn-5
	Hygromycin	500 µg/ml	Invivogen: ant-hm-5
Dissociation	Sterile PBS	-	Hyclone: SH3025601
	0.25% Trypsin-EDTA	-	Hyclone: SH3004201
	Basal Medium (see above)	40%	-
CryoMedium	Fetal Bovine Serum (FBS)	50%	Hyclone: SH3007003
	Dimethyl Sulfoxide (DMSO)	10%	Sigma: D2650

Cell handling

1. Upon receipt, directly place cells in liquid nitrogen storage. Consistent cryopreservation is essential for culture integrity.
2. Prepare Basal Medium. Prepare 37°C Water Bath. Thaw cells rapidly by removing from liquid nitrogen, and immediately immersing in a 37°C water bath, until 90% thawed. Immediately sterilize the exterior of the vial with 70% ethanol.
3. Add vial contents to 15 mL Basal Medium in T75 Tissue Culture Treated Flask. Gently swirl flask and place in a humidified, tissue culture incubator, 37°C, 5% CO₂.
4. 18-24 Hours Post–Thaw, all live cells should be attached. Viability of the cells is expected to be 60-90%, at this time, exchange Basal Medium with Selection Medium.
5. When cells are approximately 80% confluent, passage the cells. It is suggested that user expand culture to create >20 vial Master Cell Bank at low passage number. *Cells should be maintained at less than 80% confluency for optimal assay results.*
6. Cell Dissociation: Aspirate Culture Medium. Gently wash with 1x Volume PBS. Add 0.1x Volume Warm Trypsin-EDTA. Incubate 4 min, 37°C, until cells dislodge. *If cells do not round up, place in 37° C incubator for additional 2 min.* Neutralize Trypsin and collect cells in 1x Volume Basal Medium.
7. Seed Cells for expansion of culture. It is recommended that cell lines are passaged at least once before use in assays.

Table 3. Cell Culture Seeding Suggestions: *User should define based on research needs.*

Flask Size (cm ²)	Volume (mL)	Total Cell Number (x10 ⁶)	Growth Period (hrs)
T75	15	5.0	24
T75	15	2.0	48
T75	15	0.45	72

ASSAY SETUP

Fluorescence

Table 4. Settings for FLIPR^{TETRA}® with ICCD camera option

Option	Setting
Read Mode	Fluorescence
Ex/Em	Ex470_495 / Em515_575
Camera Gain	2000
Gate Open	6 %
Exposure Time	0.53
Read Interval	1s
Dispense Volume	50 µl (25 µl for 384-well)
Dispense Height	95 µl (50 µl for 384-well)
Dispense Speed	50 µl L/sec
Expel Volume	0 µl
Analysis	Subtract Bias Sample 1

Table 5. Assay Materials (Not provided)

Description	Supplier and Product Number
HBSS	Invitrogen: 14025
HEPES 1M Stock	EMD Millipore: TMS-003-C
Probenicid	Sigma:P8761
Quest Fluo-8 ^{1M} , AM	AAT Bioquest: 21080
GABA ligand	Tocris: 0344
Non-Binding 96/384 well Plates (for ligand prep)	Corning: 3605/ 3574
Black (clear Bottom) cell assay plates	Corning: 3904/ 3712

Assay Protocol - Fluorescence

1. Dissociate Culture as Recommended. Collect in Basal Medium. Document Cell Count and Viability
2. Centrifuge the cell suspension at 190 x g for six min
3. Remove supernatant. Gently resuspend the cell pellet in Basal Medium. *It is suggested that end user optimize cell plating based on individual formats.* (Default: Resuspend in volume to achieve 5×10^5 cells/ml (i.e, if collected 5×10^6 TC, $\frac{5 \times 10^6}{5 \times 10^5/ml} = 10$ mL volume)
4. Seed cell suspension into black, clear bottom plate (100 μ L/well for 96-well plate). *When seeding is complete, place the assay plate at room temperature for 30 min.*
5. Move assay plate to a humidified 37°C 5% CO₂ incubator for 18-24 h.
6. Next day, prepare Assay buffer (HBSS, 20mM HEPES, 2.5 mM Probenicid, pH 7.4) and Loading buffer (Assay buffer with 5 mM Fluo8 Dye). *Note: Please prepare Fluo8 stock according to Manufacturer's Recommendations*
7. Remove medium from assay plate and wash 1X with Assay Buffer.
8. Add Loading buffer to assay plate (100 μ L/well for 96-well plate). Incubate plate for 1.5 h at room temperature, protected from light.
9. Prepare ligands in assay buffer at 3x final concentration in non-binding plates. Use Buffer Only Control Wells for Background Subtraction.
10. Create protocol for ligand addition. Please refer to FLIPR^{TETRA}® settings provided in Table 2. Set time course for 180 s, with ligand addition at 10 s.
11. After the run is complete, apply subtract bias on sample 1. We recommend using Negative Control Correction with Buffer Only Wells. Export data to according to research needs. For most Calcium Flux analysis using Export of Max Signal to end of run is sufficient.

HOST CELL

Chem-1, an adherent cell line expressing the promiscuous G-protein, G α 15.

EXOGENOUS GENE EXPRESSION

GABBR1 & GABBR2 cDNA (Accession Number: NM_021903 & NM_005458, respectively; see CODING SEQUENCE below) expressed from a proprietary PHS plasmid.

CODING SEQUENCE
GABBR1

```

                                ATG GGG CCC GGG GCC CCT TTT GCC CGG GTG GGG TGG CCA CTG CCG
                                M   G   P   G   A   P   F   A   R   V   G   W   P   L   P

CTT CTG GTT GTG ATG GCG GCA GGG GTG GCT CCG GTG TGG GCC TCC CAC TCC CCC CAT CTC CCG CGG CCT
L   L   V   V   M   A   A   G   V   A   P   V   W   A   S   H   S   P   H   L   P   R   P

CAC TCG CGG GTC CCC CCG CAC CCC TCC TCA GAA CGG CGC GCA GTG TAC ATC GGG GCA CTG TTT CCC ATG
H   S   R   V   P   P   H   P   S   S   E   R   R   A   V   Y   I   G   A   L   F   P   M

AGC GGG GGC TGG CCA GGG GGC CAG GCC TGC CAG CCC GCG GTG GAG ATG GCG CTG GAG GAC GTG AAT AGC
S   G   G   W   P   G   G   Q   A   C   Q   P   A   V   E   M   A   L   E   D   V   N   S

CGC AGG GAC ATC CTG CCG GAC TAT GAG CTC AAG CTC ATC CAC CAC GAC AGC AAG TGT GAT CCA GGC CAA
R   R   D   I   L   P   D   Y   E   L   K   L   I   H   H   D   S   K   C   D   P   G   Q

GCC ACC AAG TAC CTA TAT GAG CTG CTC TAC AAC GAC CCT ATC AAG ATC ATC CTT ATG CCT GGC TGC AGC
A   T   K   Y   L   Y   E   L   L   Y   N   D   P   I   K   I   I   L   M   P   G   C   S

TCT GTC TCC ACG CTG GTG GCT GAG GCT GCT AGG ATG TGG AAC CTC ATT GTG CTT TCC TAT GGC TCC AGC
S   V   S   T   L   V   A   E   A   A   R   M   W   N   L   I   V   L   S   Y   G   S   S

TCA CCA GCC CTG TCA AAC CGG CAG CGT TTC CCC ACT TTC TTC CGA ACG CAC CCA TCA GCC ACA CTC CAC
S   P   A   L   S   N   R   Q   R   F   P   T   F   F   R   T   H   P   S   A   T   L   H

AAC CCT ACC CGC GTG AAA CTC TTT GAA AAG TGG GGC TGG AAG AAG ATT GCT ACC ATC CAG CAG ACC ACT
N   P   T   R   V   K   L   F   E   K   W   G   W   K   K   I   A   T   I   Q   Q   T   T

GAG GTC TTC ACT TCG ACT CTG GAC GAC CTG GAG GAA CGA GTG AAG GAG GCT GGA ATT GAG ATT ACT TTC
E   V   F   T   S   T   L   D   D   L   E   E   R   V   K   E   A   G   I   E   I   T   F

CGC CAG AGT TTC TTC TCA GAT CCA GCT GTG CCC GTC AAA AAC CTG AAG CGC CAG GAT GCC CGA ATC ATC
R   Q   S   F   F   S   D   P   A   V   P   V   K   N   L   K   R   Q   D   A   R   I   I

GTG GGA CTT TTC TAT GAG ACT GAA GCC CGG AAA GTT TTT TGT GAG GTG TAC AAG GAG CGT CTC TTT GGG
V   G   L   F   Y   E   T   E   A   R   K   V   F   C   E   V   Y   K   E   R   L   F   G

AAG AAG TAC GTC TGG TTC CTC ATT GGG TGG TAT GCT GAC AAT TGG TTC AAG ATC TAC GAC CCT TCT ATC
K   K   Y   V   W   F   L   I   G   W   Y   A   D   N   W   F   K   I   Y   D   P   S   I

AAC TGC ACA GTG GAT GAG ATG ACT GAG GCG GTG GAG GGC CAC ATC ACA ACT GAG ATT GTC ATG CTG AAT
N   C   T   V   D   E   M   T   E   A   V   E   G   H   I   T   T   E   I   V   M   L   N

CCT GCC AAT ACC CGC AGC ATT TCC AAC ATG ACA TCC CAG GAA TTT GTG GAG AAA CTA ACC AAG CGA CTG
P   A   N   T   R   S   I   S   N   M   T   S   Q   E   F   V   E   K   L   T   K   R   L

AAA AGA CAC CCT GAG GAG ACA GGA GGC TTC CAG GAG GCA CCG CTG GCC TAT GAT GCC ATC TGG GCC TTG
K   R   H   P   E   E   T   G   G   F   Q   E   A   P   L   A   Y   D   A   I   W   A   L

GCA CTG GCC CTG AAC AAG ACA TCT GGA GGA GGC GGC CGT TCC GGC GTG CGC CTG GAG GAC TTC AAC TAC
A   L   A   L   N   K   T   S   G   G   G   G   R   S   G   V   R   L   E   D   F   N   Y

AAC AAC CAG ACC ATT ACC GAC CAA ATC TAC CGG GCA ATG AAC TCC TCG TCC TTT GAG GGT GTC TCT GGC
N   N   Q   T   I   T   D   Q   I   Y   R   A   M   N   S   S   S   F   E   G   V   S   G

CAT GTG GTG TTT GAT GCC AGC GGC TCT CGG ATG GCA TGG ACG CTT ATC GAG CAG CTT CAG GGT GGC AGC
H   V   V   F   D   A   S   G   S   R   M   A   W   T   L   I   E   Q   L   Q   G   G   S

TAC AAG AAG ATT GGC TAC TAT GAC AGC ACC AAG GAT GAT CTT TCC TGG TCC AAA ACA GAT AAA TGG ATT
Y   K   K   I   G   Y   Y   D   S   T   K   D   D   L   S   W   S   K   T   D   K   W   I

GGA GGG TCC CCC CCA GCT GAC CAG ACC CTG GTC ATC AAG ACA TTC CGC TTC CTG TCA CAG AAA CTC TTT
G   G   S   P   P   A   D   Q   T   L   V   I   K   T   F   R   F   L   S   Q   K   L   F

ATC TCC GTC TCA GTT CTC TCC AGC CTG GGC ATT GTC CTA GCT GTT GTC TGT CTG TCC TTT AAC ATC TAC
I   S   V   S   V   L   S   S   L   G   I   V   L   A   V   V   C   L   S   F   N   I   Y

AAC TCA CAT GTC CGT TAT ATC CAG AAC TCA CAG CCC AAC CTG AAC AAC CTG ACT GCT GTG GGC TGC TCA
N   S   H   V   R   Y   I   Q   N   S   Q   P   N   L   N   N   L   T   A   V   G   C   S

CTG GCT TTA GCT GCT GTC TTC CCC CTG GGG CTC GAT GGT TAC CAC ATT GGG AGG AAC CAG TTC CCT TTC

```

L A L A A V F P L G L D G Y H I G R N Q F P F
 GTC TGC CAG GCC CGC CTC TGG CTC CTG GGC CTG GGC TTT AGT CTG GGC TAC GGT TCC ATG TTC ACC AAG
 V C Q A R L W L L G L G F S L G Y G S M F T K
 ATT TGG TGG GTC CAC ACG GTC TTC ACA AAG AAG GAA GAA AAG AAG GAG TGG AGG AAG ACT CTG GAA CCC
 I W W V H T V F T K K E E K K E W R K T L E P
 TGG AAG CTG TAT GCC ACA GTG GGC CTG CTG GTG GGC ATG GAT GTC CTC ACT CTC GCC ATC TGG CAG ATC
 W K L Y A T V G L L V G M D V L T L A I W Q I
 GTG GAC CCT CTG CAC CGG ACC ATT GAG ACA TTT GCC AAG GAG GAA CCT AAG GAA GAT ATT GAC GTC TCT
 V D P L H R T I E T F A K E E P K E D I D V S
 ATT CTG CCC CAG CTG GAG CAT TGC AGC TCC AGG AAG ATG AAT ACA TGG CTT GGC ATT TTC TAT GGT TAC
 I L P Q L E H C S S R K M N T W L G I F Y G Y
 AAG GGG CTG CTG CTG CTG CTG GGA ATC TTC CTT GCT TAT GAG ACC AAG AGT GTG TCC ACT GAG AAG ATC
 K G L L L L L G I F L A Y E T K S V S T E K I
 AAT GAT CAC CGG GCT GTG GGC ATG GCT ATC TAC AAT GTG GCA GTC CTG TGC CTC ATC ACT GCT CCT GTC
 N D H R A V G M A I Y N V A V L C L I T A P V
 ACC ATG ATT CTG TCC AGC CAG CAG GAT GCA GCC TTT GCC TTT GCC TCT CTT GCC ATA GTT TTC TCC TCC
 T M I L S S Q Q D A A F A F A S L A I V F S S
 TAT ATC ACT CTT GTT GTG CTC TTT GTG CCC AAG ATG CGC AGG CTG ATC ACC CGA GGG GAA TGG CAG TCG
 Y I T L V V L F V P K M R R L I T R G E W Q S
 GAG GCG CAG GAC ACC ATG AAG ACA GGG TCA TCG ACC AAC AAC AAC GAG GAG GAG AAG TCC CGG CTG TTG
 E A Q D T M K T G S S T N N N E E E K S R L L
 GAG AAG GAG AAC CGT GAA CTG GAA AAG ATC ATT GCT GAG AAA GAG GAG CGT GTC TCT GAA CTG CGC CAT
 E K E N R E L E K I I A E K E E R V S E L R H
 CAA CTC CAG TCT CGG CAG CAG CTC CGC TCC CGG CGC CAC CCA CCG ACA CCC CCA GAA CCC TCT GGG GGC
 Q L Q S R Q Q L R S R R H P P T P P E P S G G
 CTG CCC AGG GGA CCC CCT GAG CCC CCC GAC CGG CTT AGC TGT GAT GGG AGT CGA GTG CAT TTG CTT TAT
 L P R G P P E P P D R L S C D G S R V H L L Y
 AAG TGA
 K Stp

GABBR2

ATG GCT TCC CCG CCT CCC GCA CTC AGC TCG CTC CCA
 M A S P P P A L S S L P
 CCC CTT CCC GGC GTG ATT GGT CCG TCA CGG GCG CCG CCG CCG CCA CCG CCG CCC GCG CGC CTG CTA CTG
 P L P G V I G P S R A P P P P P P A R L L L
 CTA CTG CTG TAG CCG CTG CTG CCT CTG CCG CCC GGG GCC TGG GGC TGG GCG CGG GGC GCC CCC CTG
 L L L L P L L L P L A P G A W G ▼ W A R G A P R
Predicted native signal sequence protease cleavage site
 CCG CCG CCC AGC AGC CCG CCG CTC TCC ATC ATG GGC CTC ATG CCG CTC ACC AAG GAG GTG GCC AAG GGC
 P P P S S P P L S I M G L M P L T K E V A K G
 AGC ATC GGG CGC GGT GTG CTC CCC GCC GTG GAA CTG GCC ATC GAG CAG ATC CGC AAC GAG TCA CTC CTG
 S I G R G V L P A V E L A I E Q I R N E S L L
 CGC CCC TAC TTC CTC GAC CTG CGG CTC TAT GAC ACG GAG TGC GAC AAC GCA AAA GGG TTG AAA GCC TTC
 R P Y F L D L R L Y D T E C D N A K G L K A F
 TAC GAT GCA ATA AAA TAC GGG CCG AAC CAC TTG ATG GTG TTT GGA GGC GTC TGT CCA TCC GTC ACA TCC
 Y D A I K Y G P N H L M V F G V C P S V T S
 ATC ATT GCA GAG TCC CTC CAA GGC TGG AAT CTG GTG CAG CTT TCT TTT GCT GCA ACC ACG CCT GTT CTA
 I I A E S L Q G W N L V Q L S F A A T T P V L
 GCC GAT AAG AAA AAA TAC CCT TAT TTC TTT CGG ACC GTC CCA TCA GAC AAT GCG GTG AAT CCA GCC ATT
 A D K K K Y P Y F F R T V P S D N A V N P A I
 CTG AAG TTG CTC AAG CAC TAC CAG TGG AAG CGC GTG GGC ACG CTG ACG CAA GAC GTT CAG AGG TTC TCT
 L K L L K H Y Q W K R V G T L T Q D V Q R F S
 GAG GTG CGG AAT GAC CTG ACT GGA GTT CTG TAT GGC GAG GAC ATT GAG ATT TCA GAC ACC GAG AGC TTC
 E V R N D L T G V L Y G E D I E I S D T E S F

TCC AAC GAT CCC TGT ACC AGT GTC AAA AAG CTG AAG GGG AAT GAT GTG CGG ATC ATC CTT GGC CAG TTT
S N D P C T S V K K L K G N D V R I I L G Q F
GAC CAG AAT ATG GCA GCA AAA GTG TTC TGT TGT GCA TAC GAG GAG AAC ATG TAT GGT AGT AAA TAT CAG
D Q N M A A K V F C C A Y E E N M Y G S K Y Q
TGG ATC ATT CCG GGC TGG TAC GAG CCT TCT TGG TGG GAG CAG GTG CAC ACG GAA GCC GAC TCA TCC CGC
W I I P G W Y E P S W W E Q V H T E A D S S R
TGC CTC CGG AAG AAT CTG CTT GCT GCC ATG GAG GGC TAC ATT GGC GTG GAT TTC GAG CCC CTG AGC TCC
C L R K A N L L A A M E G Y I G V D F E P L S S
AAG CAG ATC AAG ACC ATC TCA GGA AAG ACT CCA CAG TAT GAG AGA GAG TAC AAC AAC AAG CGG TCA
K Q I K T I S G K T P Q Q Y E R E Y N N K R S
GGC GTG GGG CCC AGC AAG TTC CAC GGG TAC GCC TAC GAT GGC ATC TGG GTC ATC GCC AAG ACA CTG CAG
G V G P S K F H G Y A Y D G I W V I A K T L Q
AGG GCC ATG GAG ACA CTG CAT GCC AGC AGC CGG CAC CAG CGG ATC CAG GAC TTC AAC TAC ACG GAC CAC
R A M E T L H A S S R G H Q R I Q D F N Y T D H
ACG CTG GGC AGG ATC CTC AAT GSC ATG AAC GAG ACC AAC TTC TTC GGG GTC ACG GGT CAA GTT GTA
T L G R I I L N A M N E T N F F G V T G Q V V
TTC CGG AAT GGG GAG AGA ATG GGG ACC ATT AAA TTT ACT CAA TTT CAA GAC AGC AGG GAG GTG AAG GTG
F R N G E R M G T I K F T Q F Q D S R E V K V
GGA GAG TAC AAC GCT GTG GCC GAC ACA CTG GAG ATC ATC AAT GAC ACC ATC AGG TTC CAA GGA TCC GAA
G E Y N A V A D T L E I I N D T I R F Q G S E
CCA CCA AAA GAC AAC ATC ATC CTG GAG CAG CTG CGG AAG ATC TCC CTA CCT CTC TAC AGC ATC CTC
P P K D K T I I L E Q L R K I S L P L Y S I L
TCT GCC CTC ACC ATC CTC GGG ATG ATC ATG GCC AGT GCT TTT CTC TTC TTC AAC ATC AAG AAC CGG AAT
S A L T I L G M I M A S A F L F F N I K N R N
CAG AAG CTC ATA AAG ATG TCG AGT CCA TAC ATG AAC AAC CTT ATC ATC CTT GGA GGG ATG CTC TCC TAT
Q K L I K M S S P Y M N N L I I L G G M L S Y
GCT TCC ATA TTT CTC TTT GGC CTT GAT GGA TCC TTT GTC TCT GAA AAG ACC TTT GAA ACA CTT TGC ACC
A S I F L F G L D G S F V S E K T F E T L C T
GTC AGG ACC TGG ATT CTC ACC GTG GGC TAC ACG ACC GCT TTT GGG GCC ATG TTT GCA AAG ACC TGG AGA
V R T W I L T V G Y T T A F G A M F A K T W R
GTC CAC GCC ATC TTC AAA AAT GTG AAA ATG AAG AAG AAG ATC ATC AAG GAC CAG AAA CTG CTT GTG ATC
V H A I F K N V K M K K I I K D Q K L V I
GTG GGG GGC ATG CTG CTG ATC GAC CTG TGT ATC CTG ATC TGC TGG CAG GCT GTG GAC CCC CTG CGA AGG
V G G M L L I D L C I L I C W Q A V D P L R R
ACA GTG GAG AAG TAC AGC ATG GAG CCG GAC CCA GCA GGA CGG GAT ATC TCC ATC CGC CCT CTC CTG GAG
T V E K Y S M E P D P A G R D I S I R P L L E
CAC TGT GAG AAC ACC CAT ATG ACC ATC TGG CTT GGC ATC GTC TAT GCC TAC AAG GGA CTT CTC ATG TTG
H C E N T H M T I W L G I V Y A Y K G L L M L
TTC GGT TGT TTC TTA GCT TGG GAG ACC CGC AAC GTC AGC ATC CCC GCA CTC AAC GAC AGC AAG TAC ATC
F G C F L A W E T R N V S I P A L N D S K Y I
GGG ATG AGT GTC TAC AAC GTG GGG ATC ATG TGC ATC ATC GGG GCC GGT GTC TCC TTC CTG ACC CGG GAC
G M S V Y N V G I M C I I G A A V S F L T R D
CAG CCC AAT GTG CAG TTC TGC ATC GTG GCT CTG GTC ATC ATC TTC TGC AGC ACC ATC ACC CTC TGC CTG
Q P N V Q F C I V A L V I I F C S T I T L C L
GTA TTC GTG CCG AAG CTC ATC ACC CTG AGA ACA AAC CCA GAT GCA GCA ACG CAG AAC AGG CGA TTC CAG
V F V P K L I T L R T N P D A A T Q N R R F Q
TTC ACT CAG AAT CAG AAG AAA GAA GAT TCT AAA ACG TCC ACC TCG GTC ACC AGT GTG AAC CAA GCC AGC
F T Q N Q K K E D S K T S T S V T S V N Q A S
ACA TCC CGC CTG GAG GGC CTA CAG TCA GAA AAC CAT CAC CTG CGA ATG AAG ATC ACA GAG CTG GAT AAA
T S R L E G L Q S E N H H L R M K I T E L D K
GAC TTG GAA GAG GTC ACC ATG CAG CTG CAG GAC ACA CCA GAA AAG ACC ACC TAC ATT AAA CAG AAC CAC
D L E E V T M Q L Q D T P E K T T Y I K Q N H
TAC CAA GAG CTC AAT GAC ATC CTC AAC CTG GGA AAC TTC ACT GAG AGC ACA GAT GGA GGA AAG GCC ATT
Y Q E L N D I L N L G N F T E S T D G G K A A I
TTA AAA AAT CAC CTC GAT CAA AAT CCC CAG CTG CAG TGG AAC ACA ACA GAG CCC TCT CGA ACA TGC AAA
L K N H L D Q N P Q L Q W N T T E P S R T C K
GAT CCT ATA GAA GAT ATA AAC TCT CCA GAA CAC ATC CAG CGT CGG CTG TCC CTC CAG CTC CCC ATC CTC
D P I E D I N S P E H I Q R R L S L Q L P I L
CAC CAC GCC TAC CTC CCA TCC ATC GGA GGC GTG GAC GCC AGC TGT GTC AGC CCC TGC GTC AGC CCC ACC
H H A Y L P S I G G V D A S C V S P C V S P T
GCC AGC CCC CGC CAC AGA CAT GTG CCA CCC TCC TTC CGA GTC ATG GTC TCG GGC CTG TGA
A S P R H R H V P P S F R V M V S G L Stp

RELATED PRODUCTS

Product Number	Description
HTSCHEM-1	ChemiScreen™ Chem-1 parental cells
HTS119M	ChemiScreen™ GABA _B GABA Family Receptor membrane prep
HTS119RTA	Ready-to-Assay™ GABA _B GABA Family Receptor frozen cells

REFERENCES

1. Bowery NG *et al.* (2002) International Union of Pharmacology. XXXIII. Mammalian γ -aminobutyric acid_B receptors: Structure and function. *Pharmacol. Rev.* 54: 247-264.

FOR RESEARCH USE ONLY; NOT FOR USE IN DIAGNOSTIC PROCEDURES. NOT FOR HUMAN OR ANIMAL CONSUMPTION

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

No part of these works may be reproduced in any form without permission in writing.

Limited Use Label License Agreement

In addition to the General Terms & Conditions of Sale for Products and Services section, this Product is subject to Limited Use Label License Agreement. Please go to <https://www.eurofinsdiscoveryservices.com/cms/cms-content/misc/legal-disclaimer/> for more information.

Eurofins Pharma Bioanalytics Services US Inc. is an independent member of Eurofins Discovery Services