

PRODUCT DATASHEET

ChemiScreen™ CaS Calcium Sensor Receptor Stable Cell Line

CATALOG NUMBER: HTS137C

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.

Calcium homeostasis in vertebrates is regulated by hormonal communication between the parathyroid and thyroid glands, kidney, bone, and intestine. A central mediator of this regulatory system is the calcium sensor CaS (also known as CaR), a class III GPCR that is activated by high concentrations of extracellular calcium and other divalent cations. The cell types that are responsive to variations in blood calcium concentration express CaS and respond to changes in extracellular calcium in a CaS-dependent fashion. Activation of CaS in parathyroid chief cells inhibits PTH release, which results in decreased calcium mobilization from bone. In contrast, stimulation of CaS activity in the C cells of the thyroid increases secretion of calcitonin, which causes decreased bone resorption and increased urinary excretion of calcium. Allosteric "calcimimetic" potentiators are being evaluated for treatment of primary hyperparathyroidism (Brown and MacLeod, 2001). The cloned human CaS-expressing cell line is made in the Chem-1 host, which supports high levels of recombinant CaS expression on the cell surface and contains high levels of the promiscuous G protein Gα15 to couple the receptor to the calcium signaling pathway. Thus, the cell line is an ideal tool for screening for antagonists of interactions between CaS and its ligands.

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

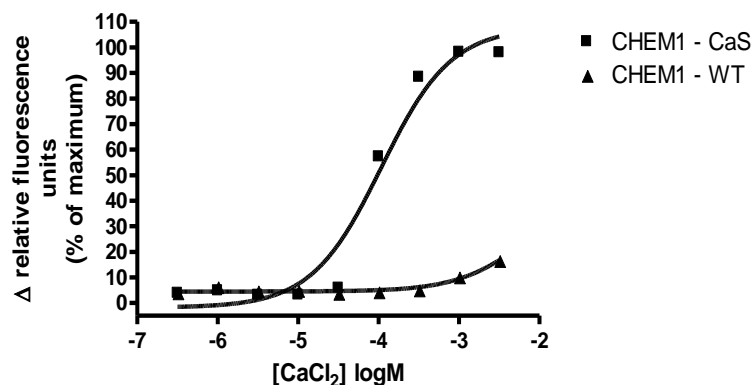


Figure 1. Representative data for activation of CaS receptor stably expressed in Chem-1 cells induced by CaCl₂ using a fluorescent calcium flux assay. CaS-expressing Chem-1 cells were seeded at 50,000 cells per well into a 96-well plate, and the following day the cells were loaded with a fluorescent 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 6,000 RLU. Similarly parental cells (catalog #: HTSCHEM-1) were tested to determine the specificity of the resulting signal.

Table 1. EC₅₀ value of CaS-expressing Chem-1 cells.

LIGAND	ASSAY	POTENCY EC ₅₀ (nM)	REFERENCE
CaCl ₂	Calcium Flux - Fluorescence	11	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.

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: SH30070.03
	Non-Essential Amino Acids (NEAA)	1X	Hyclone: SH30238.01
	HEPES	1X	EMD Millipore: TMS-003-C
Selection Medium	Basal Medium (see above)	-	
	Geneticin (G418)	250 µg/ml	Invivogen: ant-gn-5
Dissociation	Sterile PBS	-	Hyclone: SH30028.03
	0.25% Trypsin-EDTA	-	Hyclone: SH30042.01
CryoMedium	Basal Medium (see above)	40%	
	Fetal Bovine Serum (FBS)	50%	Hyclone: SH30070.03
	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/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™, AM	AAT Bioquest: 21080
CaCl ₂	Sigma: C1016
Non-Binding 96/384 well Plates (for ligand prep)	Corning: 3605/ 3574
Black (clear Bottom) cell assay plates	Corning: 3904/ 3712
Coelenterazine-h (250µg). Prepare to 10mM	Promega: S2011

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 5x10⁵ cells/ml (i.e, if collected 5e6 TC, ^{5e6/5e5/ml} = 10 mL volume)
4. Seed cell suspension into black, clear bottom plate (100 µL/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 µL/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, Gα15.

EXOGENOUS GENE EXPRESSION

Human CaS cDNA (Accession Number: NM_000388; see CODING SEQUENCE below) and promiscuous G protein are expressed in a bicistronic vector

CODING SEQUENCE

ATG GCA TTT TAT AGC TGC TGC TGG GTC CTC TTG GCA CTC ACC TGG CAC ACC TCT GCC CAC GGG CCA GAC	69
M A F Y S C C W V L L A L T W H T S A H G P D	23
CAG CGA GCC CAA AAG AAG GGG GAC ATT ATC CTT GGG GGG CTC TTT CCT ATT CAT TTT GGA GTA GCA GCT	138
Q R A Q K K G D I I L G G L F P I H F G V A A	46
AAA GAT CAA GAT CTC AAA TCA AGG CCG GAG TCT GTG GAA TGT ATC AGG TAT AAT TTC CGT GGG TTT CGC	207
K D Q D L K S R P E S V E C I R Y N F R G F R	69
TGG TTA CAG GCT ATG ATA TTT GCC ATA GAG GAG ATA AAC AGC AGC CCA GCC CTT CTT CCC AAC TTG ACG	276
W L Q A M I F A I E E I N S S P A L L P N L T	92
CTG GGA TAC AGG ATA TTT GAC ACT TGC AAC ACC GTT TCT AAG GCC TTG GAA GCC ACC CTG AGT TTT GTT	345
L G Y R I F D T C N T V S K A L E A T L S F V	115
GCT CAA AAC AAA ATT GAT TCT TTG AAC CTT GAT GAG TTC TGC AAC TGC TCA GAG CAC ATT CCC TCT ACG	414
A Q N K I D S L N L D E F C N C S E H I P S T	138
ATT GCT GTG GTG GGA GCA ACT GGC TCA GGC GTC TCC ACG GCA GTG GCA AAT CTG CTG GGG CTC TTC TAC	483
I A V V G A T G S G V S T A V A N L L G L F Y	161
ATT CCC CAG GTC AGT TAT GCC TCC TCC AGC AGA CTC CTC AGC AAC AAG AAT CAA TTC AAG TCT TTC CTC	552
I P Q V S Y A S S S R L L S N K N Q F K S F L	184
CGA ACC ATC CCC AAT GAT GAG CAC CAG GCC ACT GCC ATG GCA GAC ATC ATC GAG TAT TTC CGC TGG AAC	621
R T I P N D E H Q A T A M A D I I E Y F R W N	207
TGG GTG GGC ACA ATT GCA GCT GAT GAC GAC TAT GGG CGG CCG GGG ATT GAG AAA TTC CGA GAG GAA GCT	690
W V G T I A A D D D Y G R P G I E K F R E E A	230
GAG GAA AGG GAT ATC TGC ATC GAC TTC AGT GAA CTC ATC TCC CAG TAC TCT GAT GAG GAA GAG ATC CAG	759
E E R D I C I D F S E L I S Q Y S D E E E I Q	253
CAT GTG GTA GAG GTG ATT CAA AAT TCC ACG GCC AAA GTC ATC GTG GTT TTC TCC AGT GGC CCA GAT CTT	828
H V V E V I Q N S T A K V I V V F S S G P D L	276
GAG CCC CTC ATC AAG GAG ATT GTC CGG CGC AAT ATC ACG GGC AAG ATC TGG CTG GCC AGC GAG GCC TGG	897
E P L I K E I V R R N I T G K I W L A S E A W	299
GCC AGC TCC TCC CTG ATC GCC ATG CCT CAG TAC TTC CAC GTG GTT GGC GGC ACC ATT GGA TTC GCT CTG	966
A S S S L I A M P Q Y F H V V G G T I G F A L	322
AAG GCT GGG CAG ATC CCA GGC TTC CGG GAA TTC CTG AAG AAG GTC CAT CCC AGG AAG TCT GTC CAC AAT	1035
K A G Q I P G F R E F L K K V H P R K S V H N	345
GGT TTT GCC AAG GAG TTT TGG GAA GAA ACA TTT AAC TGC CAC CTC CAA GAA GGT GCA AAA GGA CCT TTA	1104
G F A K E F W E E T F N C H L Q E G A K G P L	368
CCT GTG GAC ACC TTT CTG AGA GGT CAC GAA GAA AGT GGC GAC AGG TTT AGC AAC AGC TCG ACA GCC TTC	1173
P V D T F L R G H E E S G D R F S N S S T A F	391
CGA CCC CTC TGT ACA GGG GAT GAG AAC ATC AGC AGT GTC GAG ACC CCT TAC ATA GAT TAC ACG CAT TTA	1242
R P L C T G D E N I S S V E T P Y I D Y T H L	414
CGG ATA TCC TAC AAT GTG TAC TTA GCA GTC TAC TCC ATT GCC CAC GCC TTG CAA GAT ATA TAT ACC TGC	1311
R I S Y N V Y L A V Y S I A H A L Q D I Y T C	437
TTA CCT GGG AGA GGG CTC TTC ACC AAT GGC TCC TGT GCA GAC ATC AAG AAA GTT GAG GCG TGG CAG GTC	1380
L P G R G L F T N G S C A D I K K V E A W Q V	460
CTG AAG CAC CTA CGG CAT CTA AAC TTT ACA AAC AAT ATG GGG GAG CAG GTG ACC TTT GAT GAG TGT GGT	1449
L K H L R H L N F T N N M G E Q V T F D E C G	483
GAC CTG GTG GGG AAC TAT TCC ATC ATC AAC TGG CAC CTC TCC CCA GAG GAT GGC TCC ATC GTG TTT AAG	1518
D L V G N Y S I I N W H L S P E D G S I V F K	506
GAA GTC GGG TAT TAC AAC GTC TAT GCC AAG AAG GGA GAA AGA CTC TTC ATC AAC GAG GAG AAA ATC CTG	1587
E V G Y Y N V Y A K K G E R L F I N E E K I L	529
TGG AGT GGG TTC TCC AGG GAG GTG CCC TTC TCC AAC TGC AGC CGA GAC TGC CTG GCA GGG ACC AGG AAA	1656
W S G F S R E V P F S N C S R D C L A G T R K	552
GGG ATC ATT GAG GGG GAG CCC ACC TGC TGC TTT GAG TGT GTG GAG TGT CCT GAT GGG GAG TAT AGT GAT	1725
G I I E G E P T C C F E C V E C P D G E Y S D	575
GAG ACA GAT GCC AGT GCC TGT AAC AAG TGC CCA GAT GAC TTC TGG TCC AAT GAG AAC CAC ACC TCC TGC	1794

E	T	D	A	S	A	C	N	K	C	P	D	D	F	W	S	N	E	N	H	T	S	C	598	
ATT	GCC	AAG	GAG	ATC	GAG	TTT	CTG	TCG	TGG	ACG	GAG	CCC	TTT	GGG	ATC	GCA	CTC	ACC	CTC	TTT	GCC	GTG	1863	
I	A	K	E	I	E	F	L	S	W	T	E	P	F	G	I	A	L	T	L	F	A	V	621	
CTG	GGC	ATT	TTC	CTG	ACA	GCC	TTT	GTG	CTG	GGT	GTG	TTT	ATC	AAG	TTC	CGC	AAC	ACA	CCC	ATT	GTC	AAG	1932	
L	G	I	F	L	T	A	F	V	L	G	V	F	I	K	F	R	N	T	P	I	V	K	644	
GCC	ACC	AAC	CGA	GAG	CTC	TCC	TAC	CTC	CTC	CTC	TTC	TCC	CTG	CTC	TGC	TGC	TTC	TCC	AGC	TCC	CTG	TTC	2001	
A	T	N	R	E	L	S	Y	L	L	L	F	S	L	L	C	C	F	S	S	S	S	L	F	667
TTC	ATC	GGG	GAG	CCC	CAG	GAC	TGG	ACG	TGC	CGC	CTG	CGC	CAG	CCG	GCC	TTT	GGC	ATC	AGC	TTC	GTG	CTC	2070	
F	I	G	E	P	Q	D	W	T	C	R	L	R	Q	P	A	F	G	I	S	F	V	L	690	
TGC	ATC	TCA	TGC	ATC	CTG	GTG	AAA	ACC	AAC	CGT	GTC	CTC	CTG	GTG	TTT	GAG	GCC	AAG	ATC	CCC	ACC	AGC	2139	
C	I	S	C	I	L	V	K	T	N	R	V	L	L	V	F	E	A	K	I	P	T	S	713	
TTC	CAC	CGC	AAG	TGG	TGG	GGG	CTC	AAC	CTG	CAG	TTC	CTG	CTG	GTT	TTC	CTC	TGC	ACC	TTC	ATG	CAG	ATT	2208	
F	H	R	K	W	W	G	L	N	L	Q	F	L	L	V	F	L	C	T	F	M	Q	I	736	
GTC	ATC	TGT	GTG	ATC	TGG	CTC	TAC	ACC	GCG	CCC	CCC	TCA	AGC	TAC	CGC	AAC	CAG	GAG	CTG	GAG	GAT	GAG	2277	
V	I	C	V	I	W	L	Y	T	A	P	P	S	S	Y	R	N	Q	E	L	E	D	E	759	
ATC	ATC	TTC	ATC	ACG	TGC	CAC	GAG	GGC	TCC	CTC	ATG	GCC	CTG	GGC	TTC	CTG	ATC	GGC	TAC	ACC	TGC	CTG	2346	
I	I	F	I	T	C	H	E	G	S	L	M	A	L	G	F	L	I	G	Y	T	C	L	782	
CTG	GCT	GCC	ATC	TGC	TTC	TTT	GCC	TTC	AAG	TCC	CGG	AAG	CTG	CCG	GAG	AAC	TTC	AAT	GAA	GCC	AAG	2415		
L	A	A	I	C	F	F	F	A	F	K	S	R	K	L	P	E	N	F	N	E	A	K	805	
TTC	ATC	ACC	TTC	AGC	ATG	CTC	ATC	TTC	TTC	ATC	GTC	TGG	ATC	TCC	TTC	ATT	CCA	GCC	TAT	GCC	AGC	ACC	2484	
F	I	T	F	S	M	L	I	F	F	I	V	W	I	S	F	I	P	A	Y	A	S	T	828	
TAT	GGC	AAG	TTT	GTC	TCT	GCC	GTA	GAG	GTG	ATT	GCC	ATC	CTG	GCA	GCC	AGC	TTT	GGC	TTG	CTG	GCG	TGC	2553	
Y	G	K	F	V	S	A	V	E	V	I	A	I	L	A	A	S	F	G	L	L	A	C	851	
ATC	TTC	TTC	AAC	AAG	ATC	TAC	ATC	ATT	CTC	TTC	AAG	CCA	TCC	CGC	AAC	ACC	ATC	GAG	GAG	GTG	CGT	TGC	2622	
I	F	F	N	K	I	Y	I	I	L	F	K	P	S	R	N	T	I	E	E	V	R	C	874	
AGC	ACC	GCA	GCT	CAC	GCT	TTC	AAG	GTG	GCT	GCC	CGG	GCC	ACG	CTG	CGC	CGC	AGC	AAC	GTC	TCC	CGC	AAG	2691	
S	T	A	A	H	A	F	K	V	A	A	R	A	T	L	R	R	S	N	V	S	R	K	897	
CGG	TCC	AGC	AGC	CTT	GGA	GGC	TCC	ACG	GGA	TCC	ACC	CCC	TCC	TCC	TCC	ATC	AGC	AGC	AAG	AGC	AAC	AGC	2760	
R	S	S	S	L	G	G	S	T	G	S	T	P	S	S	S	I	S	S	K	S	N	S	920	
GAA	GAC	CCA	TTC	CCA	CAG	CCC	GAG	AGG	CAG	AAG	CAG	CAG	CAG	CCG	CTG	GCC	CTA	ACC	CAG	CAA	GAG	CAG	2829	
E	D	P	F	P	Q	P	E	R	Q	K	Q	Q	Q	P	L	A	L	T	Q	Q	E	Q	943	
CAG	CAG	CAG	CCC	CTG	ACC	CTC	CCA	CAG	CAG	CAA	CGA	TCT	CAG	CAG	CAG	CCC	AGA	TGC	AAG	CAG	AAG	GTC	2898	
Q	Q	Q	P	L	T	L	P	Q	Q	Q	R	S	Q	Q	Q	P	R	C	K	Q	K	V	966	
ATC	TTT	GGC	AGC	GGC	ACG	GTC	ACC	TTC	TCA	CTG	AGC	TTT	GAT	GAG	CCT	CAG	AAG	AAC	GCC	ATG	GCC	CAC	2967	
I	F	G	S	G	T	V	T	F	S	L	S	F	D	E	P	Q	K	N	A	M	A	H	989	
AGG	AAT	TCT	ACG	CAC	CAG	AAC	TCC	CTG	GAG	GCC	CAG	AAA	AGC	AGC	GAT	ACG	CTG	ACC	CGA	CAC	CAG	CCA	3036	
R	N	S	T	H	Q	N	S	L	E	A	Q	K	S	S	D	T	L	T	R	H	Q	P	1012	
TTA	CTC	CCG	CTG	CAG	TGC	GGG	GAA	ACG	GAC	TTA	GAT	CTG	ACC	GTC	CAG	GAA	ACA	GGT	CTG	CAA	GGA	CCT	3105	
L	L	P	L	Q	C	G	E	T	D	L	D	L	T	V	Q	E	T	G	L	Q	G	P	1035	
GTG	GGT	GGA	GAC	CAG	CGG	CCA	GAG	GTG	GAG	GAC	CCT	GAA	GAG	TTG	TCC	CCA	GCA	CTT	GTA	GTG	TCC	AGT	3174	
V	G	G	D	Q	R	P	E	V	E	D	P	E	E	L	S	P	A	L	V	V	S	S	1058	
TCA	CAG	AGC	TTT	GTC	ATC	AGT	GGT	GGA	GGC	AGC	ACT	GTT	ACA	GAA	AAC	GTA	GTG	AAT	TCA	TAA				
S	Q	S	F	V	I	S	G	G	G	S	T	V	T	E	N	V	V	N	S	Stp				

RELATED PRODUCTS

Product Number	Description
HTSCHEM-1	ChemiScreen™ Chem-1 Parental Cell Line (control cells)
HTS137M	ChemiScreen™ CaS Calcium Sensor Receptor Membrane Prep

REFERENCES

1. Brown EM and MacLeod RJ (2001) Extracellular Calcium Sensing and Extracellular Calcium Signaling. *Physiol. Rev.* 81: 239-297.

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