

γ -Aminobutyric acid

γ -Aminobutyric acid (GABA) is the chief inhibitory neurotransmitter of the CNS. Reduces neuronal excitability, regulates muscle tone, relieves anxiety, promotes lean muscle growth, burns fat, stabilizes blood pressure, and improves mood. Used to reduce symptoms of premenstrual syndrome (PMS), and in treating attention deficit hyperactivity disorder (ADHD). Agonist of the γ -aminobutyric acid receptors GABAA and GABAB.

Materials Provided

Catalog Number:	92-1218
Quantity Supplied:	1 vial(s), 10 mg (re-packaged) per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	103.12 g/mol
Source:	Synthetic
Purity:	$\geq 99\%$ by TLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	RT Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	100 mM
Reconstitution Volume:	0.970 mL
Reconstitution Solvent:	H ₂ O

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com