

L-Glutamic Acid

L-glutamic acid (glutamic acid) is a naturally produced proteinogenic non-essential amino acid that plays a role in neural activation. It is the most common excitatory neurotransmitter in the CNS. Agonist of the glutamate receptors (metabotropic, kainate, NMDA, and AMPA). May be dysregulated in nervous system disorders, such as epilepsy.

Materials Provided

Catalog Number:	92-1197
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:	147.13 g/mol
Source:	Synthetic
Purity:	≥ 95%
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:	0.679 mL
Reconstitution Solvent:	1N NaOH

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