

L-Theanine

Glutamate receptor ligand

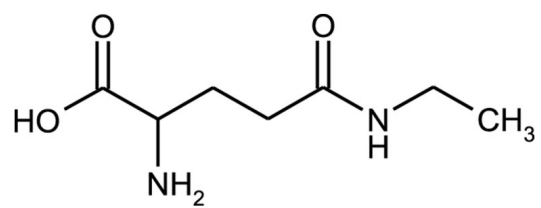
Amino acid analog of glutamine, uniquely found in green tea. Increases brain serotonin, dopamine and GABA levels and binds to AMPA, kainate and NMDA receptors. Has been shown to exert neuroprotective effects in animal models possibly through its antagonistic effects on group 1 metabotropic glutamate receptors. Behavioural studies with rats suggest L-Theanine to positively affect learning and memory processes. Promotes self-renewal of human embryonic stem cells.

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from moisture.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N γ -Ethyl-L-glutamine, L-Glutamic acid γ -(ethylamide)
Appearance	White to off-white solid.
CAS	3081-61-6
Couple Target	AMPA receptor, Kainate receptor, Metabotropic glutamate receptor, NMDA receptor
Couple Type	Inhibitor, Ligand
Formula	C ₇ H ₁₄ N ₂ O ₃
MW	174.20
Purity	≥98% (Titration)
Solubility	Soluble in water (50mg/ml).
Source	Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com