Ibotenic acid

mGluR agonist

Potent glutamate receptor agonist. Displays full agonist activity at mGlu1 and mGlu5.

Citations: 3

View Online »

Ordering Information

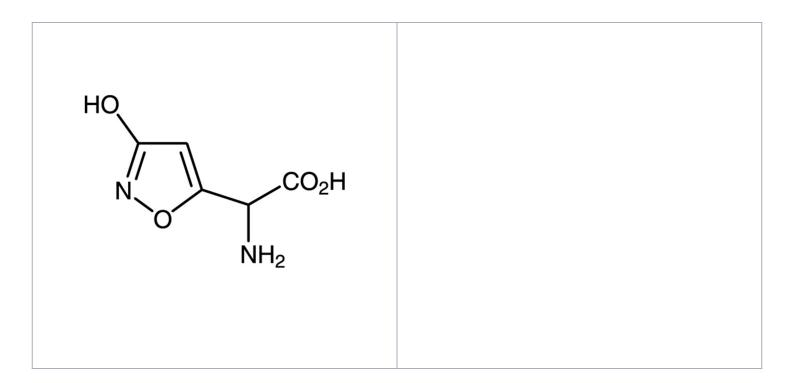
Order Online »

BML-EA120-0001

1mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied,

at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name α-Amino-3-hydroxy-5-isoxazoleacetic acid

Appearance White to off-white solid.

CAS 2552-55-8

Couple Target Metabotropic glutamate receptor

Couple Type Ligand

Formula $C_5H_6N_2O_4$

Identity Determined by 1H-NMR and MS.

MW 158.1

Purity ≥98% (HPLC)

Solublity Soluble in water (20mg/ml).

Last modified: May 29, 2024

