

Ibotenic acid

mGluR agonist

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

Citations: 3

[View Online »](#)

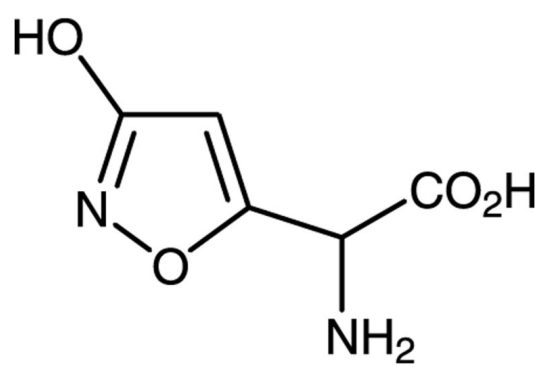
Ordering Information

[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) |
| Solubility | Soluble in water (20mg/ml). |



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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