

HOE-140

Bradykinin receptor ligand

A potent and selective bradykinin B2 antagonist. HOE-140 is a useful tool for discrimination of the physiological roles of bradykinin B2 versus B1 receptors. Displays a protective effect in myocardial reperfusion injury.

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Icatibant

Appearance White solid.

CAS 130308-48-4

Couple Target Bradykinin receptor

Couple Type Ligand

Formula $C_{59}H_{89}N_{19}O_{13}S$

Identity Determined by MS.

MW 1304.5

Purity ≥95% (HPLC)

Sequence D-Arg-Arg-Pro-Hyp-Gly-Thi-Ser-Tic-Oic-Arg

Solubility Soluble in water.

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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