

SB 225002

Chemokine receptor ligand

Potent and selective antagonist of CXCR2 chemokine receptor (IC_{50} =22nM). >150-fold more selective over CXCR1 receptor. Inhibits IL-8- (IC_{50} =8nM) and GRO α -mediated (IC_{50} =10nM) calcium mobilisation. Prevents IL-8-induced neutrophil migration. Inhibits HIV replication in lymphocytes and macrophages.

Citations: 4

[View Online »](#)

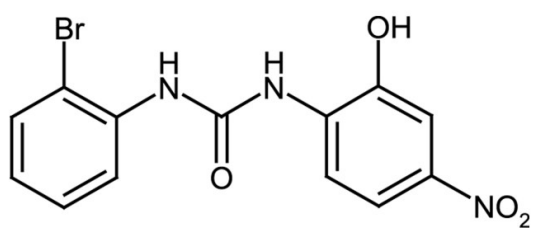
Ordering Information

[Order Online »](#)

ALX-270-478-M001	1mg
ALX-270-478-M005	5mg

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-(2-Bromophenyl)-N'-(2-hydroxy-4-nitrophenyl)urea

Appearance Yellow solid.

CAS 182498-32-4

Couple Target Chemokine receptor

Couple Type Ligand

Formula $C_{13}H_{10}BrN_3O_4$

Identity Identity determined by MS, 1H-NMR and 13C-NMR.

MW 352.1

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (40mg/ml), 100% ethanol or methanol (15mg/ml); slightly soluble in acetonitrile (2mg/ml).

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-
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)