

RHC-80267

Diacylglycerol lipase inhibitor

Diacylglycerol (DAG) is an important cellular source of arachidonate which may be released by the action of DAG lipase. RHC-80267 selectively inhibits DAG lipase activity in canine platelets ($IC_{50}=4\text{ }\mu\text{M}$), bovine adrenal chromaffin cells, human adrenal glomerulosa cells, rat thyroid lobes and pancreatic minilobules.

Citations: 8

[View Online »](#)

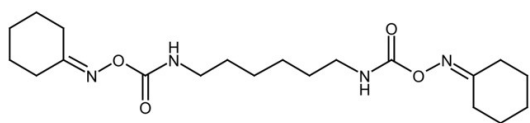
Ordering Information

[Order Online »](#)

BML-ST300-0010	10mg
BML-ST300-0050	50mg

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 2 years. 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	U-57908, 1,6-bis-(Cyclohexyloximinocarbonylamino)-hexane
Appearance	White to off-white solid.
CAS	83654-05-1
Couple Target	Diacylglycerol lipase
Couple Type	Inhibitor
Formula	$C_{20}H_{34}N_4O_4$
MW	394.6
MeltingPoint	118-121°C
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/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)