

ITSA-1

HDAC inhibitor

ITSA-1 is a cell permeable benzotriazole amide that counteracts trichostatin A (Prod. No. BML-GR309) induced cell cycle arrest, histone acetylation and transcriptional activation but not the activity of other HDAC inhibitors.

Citations: 2

[View Online »](#)

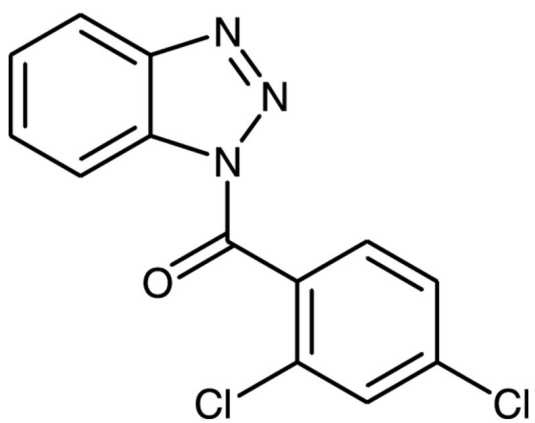
Ordering Information

[Order Online »](#)

BML-GR350-0025	25mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	N-(1H-Benzotriazol-1-yl)-2,4-dichlorobenzamide
Appearance	White solid.
CAS	200626-61-5
Couple Target	HDAC
Couple Type	Inhibitor
Formula	$C_{13}H_7Cl_2N_3O$
MW	292.1
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/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)