

BML-282

UCH inhibitor

Novel inhibitor of ubiquitin C-terminal hydrolase L1 ($IC_{50} = 0.81 \mu M$). It displays 21-fold greater selectivity over UCH-L3 ($IC_{50} = 17 \mu M$).

Citations: 1

[View Online »](#)

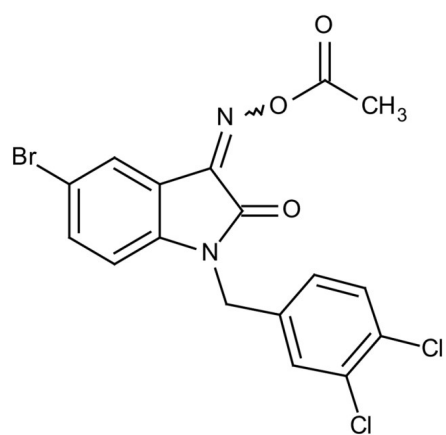
Ordering Information

[Order Online »](#)

BML-EI400-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 1 month.
Handling	Store tightly sealed. Allow product to equilibrate to room temperature prior to opening.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(3,4-Dichlorobenzyl)-5-bromoisatin O-acetyloxime
Appearance	Yellow solid.
Couple Target	Deconjugating enzyme [DCE], UCH
Couple Type	Inhibitor
Formula	$C_{17}H_{11}BrCl_2N_2O_3$
MW	442.1
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (5mg/ml with warming) or dimethyl formamide (10mg/ml with warming).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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