

TCID

UCH-L3 inhibitor

Novel inhibitor of ubiquitin C-terminal hydrolase L3 ($IC_{50} = 0.6\mu M$).
Displays 125-fold greater selectivity over UCH-L1 ($IC_{50} = 75\mu M$).

Citations: 1

[View Online »](#)

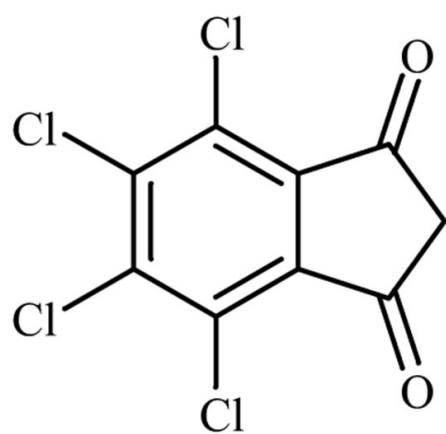
Ordering Information

[Order Online »](#)

BML-EI399-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. 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	4,5,6,7-Tetrachloroindan-1,3-dione
Appearance	Orange solid.
CAS	30675-13-9
Couple Target	Deconjugating enzyme [DCE], UCH
Couple Type	Inhibitor
Formula	$C_9H_2Cl_4O_2$
MW	283.9
Purity	≥95% (TLC)
Solubility	Soluble in DMSO (25mg/ml).

Last modified: May 29, 2024



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