

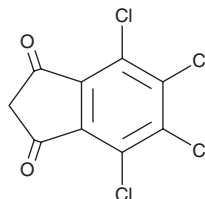
PRODUCT INFORMATION



TCID

Item No. 16353

CAS Registry No.: 30675-13-9
Formal Name: 4,5,6,7-tetrachloro-1H-indene-1,3(2H)-dione
Synonym: UCH-L3 Inhibitor
MF: $C_9H_2Cl_4O_2$
FW: 283.9
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 244 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

TCID is supplied as a crystalline solid. A stock solution may be made by dissolving the TCID in the solvent of choice, which should be purged with an inert gas. TCID is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of TCID in these solvents is approximately 10 and 15 mg/ml, respectively.

Description

Ubiquitin carboxyl-terminal hydrolase isozyme L3 (UCH-L3) is a deubiquitinating enzyme and thiol-dependent hydrolase.¹⁻⁴ TCID is a tetrachloro indandione that acts as a potent inhibitor of UCH-L3 and demonstrates 125-fold selectivity over UCH-L1 ($IC_{50}s = 0.6$ and $75 \mu\text{M}$, respectively).⁵

References

1. Mason, D.E., Ek, J., Peters, E.C., *et al.* Substrate profiling of deubiquitin hydrolases with a positional scanning library and mass spectrometry. *Biochemistry* **43**(21), 6535-6544 (2004).
2. Setsuie, R., Sakurai, M., Sakaguchi, Y., *et al.* Ubiquitin dimers control the hydrolase activity of UCH-L3. *Neurochem. Int.* **54**(5-6), 314-321 (2009).
3. Wada, H., Kito, K., Caskey, L.S., *et al.* Cleavage of the C-terminus of NEDD8 by UCH-L3. *Biochem. Biophys. Res. Commun.* **251**(3), 668-692 (1998).
4. Iphöfer, A., Kummer, A., Nimtz, M., *et al.* Profiling ubiquitin linkage specificities of deubiquitinating enzymes with branched ubiquitin isopeptide probes. *Chembiochem* **13**(10), 1416-1420 (2012).
5. Liu, Y., Lashuel, H.A., Choi, S., *et al.* Discovery of inhibitors that elucidate the role of UCH-L1 activity in the H1299 lung cancer cell line. *Chem. Biol.* **10**(9), 837-846 (2003).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/21/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM