

ML-7 . HCl

Protein kinase inhibitor

Potent and selective inhibitor of MLC kinase ($K_i=0.3 \mu\text{M}$). It inhibits also smooth-muscle myosin light chain kinase, protein kinase C (PKC) and cAMP-dependent protein kinase (PKA).

Citations: 11

[View Online »](#)

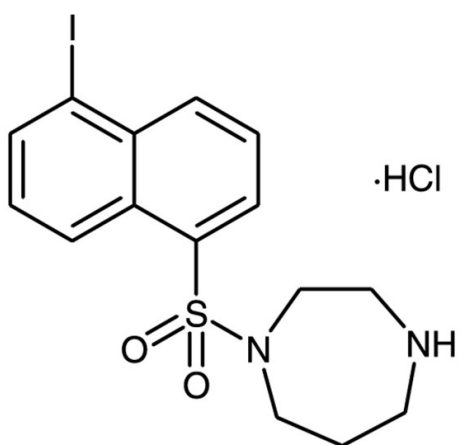
Ordering Information

[Order Online »](#)

BML-EI197-0010	10mg
----------------	------

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 -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-(5-Iodonaphthalene-1-sulfonyl)-1H-hexahydro-1,4-diazepine . HCl
Appearance	White solid.
CAS	110448-33-4
Couple Target	MLCK, PKA, PKC
Couple Type	Inhibitor
Formula	$C_{15}H_{17}IN_2O_2S \cdot HCl$
MW	452.7
Purity	≥99%
Solubility	Soluble in DMSO (10mg/ml), ethanol/water (1:1, 10mg/ml), 100% ethanol (warm) and methanol.

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