

ML-9

Myelin basic protein inhibitor

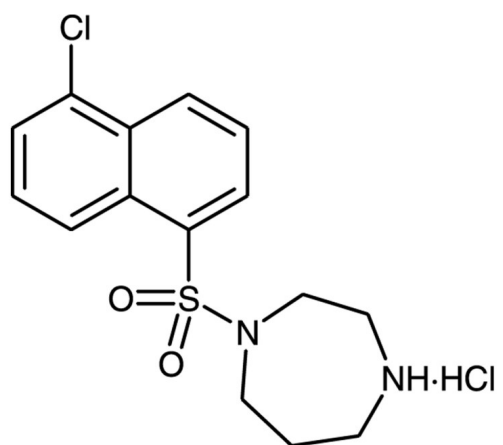
Selective inhibitor of MLC kinase ($K_i=3.8 \mu\text{M}$). Inhibits insulin-stimulated 2-deoxyglucose transport and PP-1 activation in adipocytes, possibly via an effect on MAP kinase or its activation. Useful PKB inhibitor ($\text{IC}_{50}=10\text{-}50 \mu\text{M}$). Also known to inhibit PKA ($\text{IC}_{50}=20 \mu\text{M}$) and S6 kinase ($\text{IC}_{50}=50 \mu\text{M}$). Does not inhibit PI 3-kinase.

Citations: 9

[View Online »](#)

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 0-4°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	1-(5-Chloronaphthalene-1-sulfonyl)-1H-hexahydro-1,4-diazepine . HCl
Appearance	White solid.
CAS	105637-50-1
Couple Target	Akt, MLCK, PKA, S6 kinase
Couple Type	Inhibitor
Formula	$C_{15}H_{17}N_2O_2S$ Cl . HCl
Identity	Determined by 1H-NMR and MS.
MW	361.3
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (10mg/ml) and ethanol/water (1:1, 10mg/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-be@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