

# ML-9

## Myelin basic protein inhibitor

Selective inhibitor of MLC kinase ( $K_i=3.8\text{ }\mu\text{M}$ ). 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 ( $\text{IC}_{50}=10\text{-}50\text{ }\mu\text{M}$ ). Also known to inhibit PKA ( $\text{IC}_{50}=20\text{ }\mu\text{M}$ ) and S6 kinase ( $\text{IC}_{50}=50\text{ }\mu\text{M}$ ). Does not inhibit PI 3-kinase.

Citations: 9

[View Online »](#)

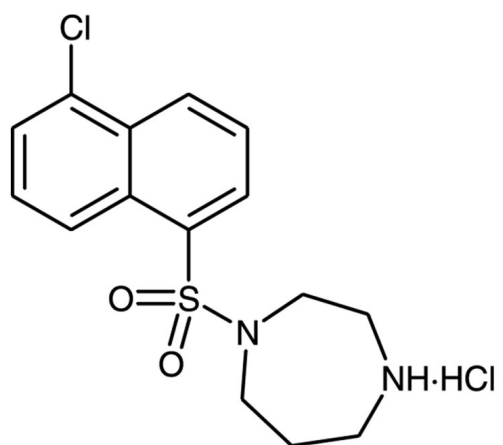
## Ordering Information

[Order Online »](#)

BML-EI153-0050	50mg
----------------	------

## 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 <sub>15</sub> H <sub>17</sub> N <sub>2</sub> O <sub>2</sub> SCI . 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](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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