

Inhibitor of trypsin and trypsin-like proteases

Irreversible inhibitor of trypsin and trypsin-like serine proteases such as granzyme D and tryptase. Also inhibits other proteases including plasmin, thrombin, and papain and some kinases including PKC. Inhibits LPS-induced expression of COX-2 in macrophages. Inhibits apoptosis (possibly through the inhibition of serine proteases). Chymotrypsin is not inhibited by TLCK, so this inhibitor can be used to block contaminating trypsin activity in chymotrypsin preparations. TLCK can be used in combination with a chymotrypsin inhibitor to prevent proteolytic degradation during the isolation of proteins.

Citations: 12

[View Online »](#)

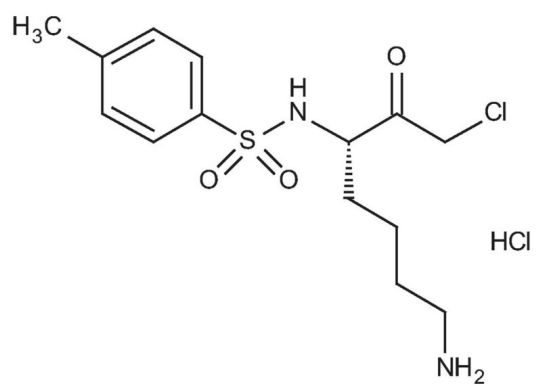
Ordering Information

[Order Online »](#)

BML-PI121-0200	200mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Na-Tosyl-Lys-chloromethylketone . HCl

Appearance Off-white solid

CAS 4238-41-9

Formula $C_{14}H_{21}ClN_2O_3S \cdot HCl$

MW 369.3

Purity ≥95% (HPLC)

Sequence Na-Tosyl-Lys-chloromethylketone

Solubility Soluble in methanol (50mg/ml), DMSO (5mM) or water. Stock solutions at concentrations up to 10mM can be prepared in aqueous buffers with a pH of 6.0 or lower. TLCK is very unstable in buffers with a pH higher than 6 and will decompose rapidly.



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)