TLCK

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 chymoptrypsin 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

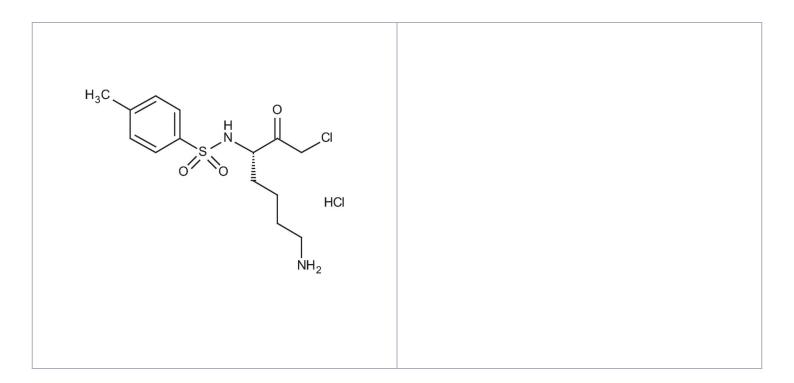
Order Online »

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 $N\alpha$ -Tosyl-Lys-chloromethylketone . HCl

Appearance Off-white solid

CAS 4238-41-9

Formula $C_{14}H_{21}CIN_2O_3S$. HCI

MW 369.3

Purity ≥95% (HPLC)

Sequence Na-Tosyl-Lys-chloromethylketone

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.

Last modified: May 29, 2024

