

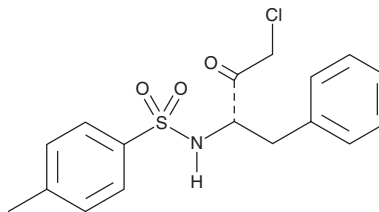
PRODUCT INFORMATION



Tosyl Phenylalanyl Chloromethyl Ketone

Item No. 20166

CAS Registry No.: 402-71-1
Formal Name: N-[(1S)-3-chloro-2-oxo-1-(phenylmethyl)propyl]-4-methyl-benzenesulfonamide
Synonyms: N^α-Tosyl-L-Phenylalanine Chloromethyl Ketone, TPCK
MF: C₁₇H₁₈ClNO₃S
FW: 351.8
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Tosyl Phenylalanyl Chloromethyl Ketone (TPCK) is supplied as a crystalline solid. A stock solution may be made by dissolving the TPCK in the solvent of choice, which should be purged with an inert gas. TPCK is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TPCK in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

TPCK is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TPCK should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. TPCK has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TPCK is an irreversible inhibitor of chymotrypsin-like proteases that has been shown to affect cell proliferation, apoptosis, and tumorigenesis.¹ It can disrupt PDK1 signaling to the AGC kinases, Akt, S6K1, and RSK, as well as MSK1 and MSK2.² TPCK also inhibits superoxide production and suppresses neutrophil respiratory burst.³

References

1. Blair, T.T. and Marini, M.A. A reexamination of the reaction of TPCK with α-chymotrypsin. *FEBS Lett.* **20(1)**, 41-43 (1972).
2. Anjum, R., Pae, E., Blenis, J., *et al.* TPCK inhibits AGC kinases by direct activation loop adduction at phenylalanine-directed cysteine residues. *FEBS Lett.* **586(19)**, 3471-3476 (2012).
3. Gillibert, M., Dehry, Z., Terrier, M., *et al.* Another biological effect of tosylphenylalanylchloromethane (TPCK): It prevents p47phox phosphorylation and translocation upon neutrophil stimulation. *Biochem J.* **386(Pt 3)** 549-556 (2005).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/26/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM