

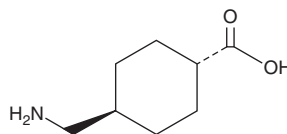
PRODUCT INFORMATION



Tranexamic Acid

Item No. 19193

CAS Registry No.: 1197-18-8
Formal Name: *trans*-4-(aminomethyl)-cyclohexanecarboxylic acid
Synonyms: AMCA, TXA
MF: C₈H₁₅NO₂
FW: 157.2
Purity: ≥95%
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥4 years



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

Laboratory Procedures

Tranexamic acid (TXA) is supplied as a crystalline solid. TXA is sparingly soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of TXA be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of TXA in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Tranexamic acid is an inhibitor of fibrinolysis that inhibits the interaction of plasmin with fibrin (IC₅₀ = 3.1 μM in isolated human plasma).¹ It decreases α-naphthylisothiocyanate (ANIT)-induced increases in the expression of the profibrogenic genes *Timp1* and *Col1a1*, but not *Itgb6*, *Tgfb1*, or *Tgfb2*, and reduces protein levels of cytokeratin 19 (CK19) and type I collagen in the liver after 28 days in a mouse model of chronic bile duct injury.²

References

1. Boström, J., Grant, J.A., Fjellström, O., *et al.* Potent fibrinolysis inhibitor discovered by shape and electrostatic complementarity to the drug tranexamic acid. *J. Med. Chem.* **56**(8), 3273-3280 (2013).
2. Joshi, N., Kopec, A.K., Towery, K., *et al.* The antifibrinolytic drug tranexamic acid reduces liver injury and fibrosis in a mouse model of chronic bile duct injury. *J. Pharmacol. Exp. Ther.* **349**(3), 383-392 (2014).

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, 12/19/2022

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