

PRODUCT INFORMATION



Tianeptine (sodium salt)

Item No. 25950

CAS Registry No.: 30123-17-2

Formal Name: 7-[(3-chloro-6,11-dihydro-6-methyl-5,5-dioxidodibenzo[c,f][1,2]thiazepin-11-yl) amino]-heptanoic acid, monosodium salt

MF: $C_{21}H_{24}ClN_2O_4S \cdot Na$

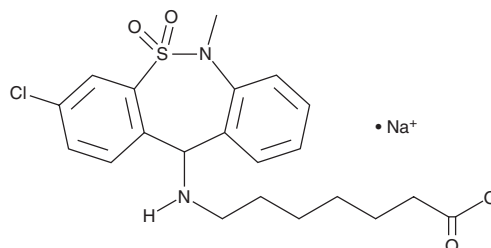
FW: 458.9

Purity: $\geq 98\%$

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 5 years



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

Description

Tianeptine (sodium salt) (Item No. 25950) is an analytical reference standard categorized as an opioid.¹ Tianeptine has been abused and associated with overdose and death.² Formulations containing tianeptine have been used in the treatment of depression.³ This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 17561).

References

1. Gassaway, M.M., Rives, M.L., Kruegel, A.C., *et al.* The atypical antidepressant and neurorestorative agent tianeptine is a μ -opioid receptor agonist. *Transl. Psychiatry* **4**, e411 (2014).
2. Lauhan, R., Hsu, A., Alam, A., *et al.* Tianeptine abuse and dependence: Case report and literature review. *Psychosomatics* **59**(6), 547-553 (2018).
3. McEwen, B.S., Chattarji, S., Diamond, D.M., *et al.* The neurobiological properties of tianeptine (Stablon): From monoamine hypothesis to glutamatergic modulation. *Mol. Psychiatry* **15**(3), 237-249 (2010).

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, 07/01/2023

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