

PRODUCT INFORMATION

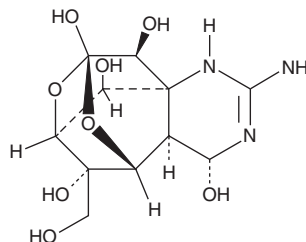


Tetrodotoxin (citrate)

Item No. 14964

CAS Registry No.: 18660-81-6
Formal Name: (4R,4aR,5R,7S,9S,10S,10aR,11S,12S)-2-amino-1,4,4a,5,9,10-hexahydro-12-(hydroxymethyl)-5,9:7,10a-dimethanol-10aH-[1,3]dioxocino[6,5-d]pyrimidine-4,7,10,11,12-pentol, citrate

Synonym: TTX citrate
MF: C₁₁H₁₇N₃O₈
FW: 319.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Tetrodotoxin (TTX) (citrate) is supplied as a solid. For biological experiments, we suggest that organic solvent-free aqueous solutions of TTX (citrate) be prepared by directly dissolving the solid in aqueous buffers. The solubility of TTX (citrate) in citrate buffer, pH 4.8, is approximately 30 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

TTX is a potent marine-derived neurotoxin that reversibly inhibits the inward sodium current through voltage-activated sodium (Na_v) channels, blocking nerve and muscle action potentials. It inhibits the Na_v current in frog muscle and squid axon with IC₅₀ values of 4.1 and 5.2 nM, respectively, and binds to rat brain membranes with a K_d value of 1.8 nM.¹ TTX was used in Hodgkin and Huxley's classic experiments to elucidate the physical biology of nerve action potentials and remains an indispensable neuroscience tool to pharmacologically dissect the contribution of the Na_v current in excitable neurons and to rationally design compounds for the treatment of neuropathic pain.¹ This citrate-stabilized form of TTX is designed for improved solubility over pure TTX (Item No. 14963), which is insoluble in water and most organic solvents.

Reference

1. Moczydlowski, E.G. The molecular mystique of tetrodotoxin. *Toxicon* **63**, 165-183 (2013).

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, 10/11/2017

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