

Tetrodotoxin

Na⁺ channel blocker

Tetrodotoxin is a non-peptide toxin isolated from ovaries and liver of Chinese puffer fish. Neurotoxin. Potent and reversible inhibitor of the voltage dependent Na⁺ channels in nerve membranes.

May require a license for import, please [contact us](#) for more information.

Citations: 27

[View Online »](#)

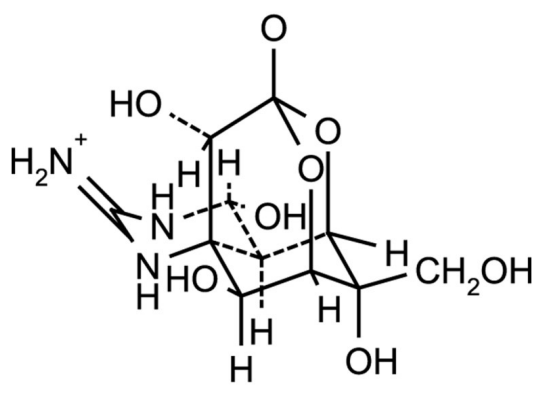
Ordering Information

[Order Online »](#)

BML-NA120-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Avoid strongly acidic or alkaline solutions. USE EXTREME CAUTION! MAY CAUSE DEATH.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TTX, Fugu poison
Appearance	White to off-white solid.
CAS	4368-28-9
Couple Target	Sodium channel
Couple Type	Blocker
Formula	$C_{11}H_{17}N_3O_8$
MI	14: 9246
MW	319.3
Purity	≥98% (HPLC)
RTECS	IO1450000

Solubility Soluble in dilute acetic acid (e.g. 1mg/ml in 10mM HOAc or 10mg/ml in 0.1M HOAc) or in citric acid/sodium citrate buffer, pH 4.8 (e.g. 1mg/ml in 20mM buffer). Do not use alkaline solution.



ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 595
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 528
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com