

# Tetrodotoxin

Na<sup>+</sup> 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<sup>+</sup> channels in nerve membranes.

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

Citations: 27

[View Online »](#)

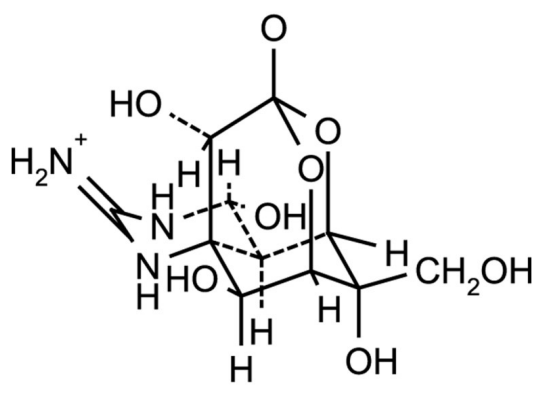
## Ordering Information

[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 <sub>11</sub> H <sub>17</sub> N <sub>3</sub> O <sub>8</sub>
MI	<b>14:</b> 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](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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