

Tetrodotoxin

This toxin reversibly blocks the sodium channels of neurons and has been a widely used tool for mechanistic studies of neurons, skeletal muscle, and cardiac muscle cells.



Product attributes

Format	Citrate Free, With Citrate
Toxin	Tetrodotoxin
CAS number	4368-28-9
Molecular weight	319.28

Product Description

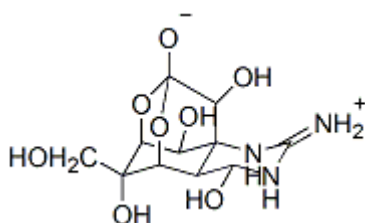
Please Note: Due to federal and state regulations, this item is can be ordered by quote request only.

Tetrodotoxin (TTX) is potent neurotoxin that reversibly binds and blocks sodium channels to disrupt the action potential and inhibit the firing of neurons. It is a widely used tool to study the mechanism of action of neurons, skeletal muscle cells, and cardiac muscle cells.

- Reversibly binds and blocks voltage-gated sodium channels
- Available with citrate or citrate-free
- Tetrodotoxin with citrate is water soluble

Tetrodotoxin citrate-free can be dissolved in slightly acidic buffer. Tetrodotoxin with citrate can be dissolved in water to obtain TTX in pH 4.8 citrate buffer.

Chemical Structure:



Note: This product is not available for sale outside of the U.S.

References

1. Neurocomputing. 70, 10 (2007), DOI: [10.1016/j.neucom.2006.10.135](https://doi.org/10.1016/j.neucom.2006.10.135)
2. Frontiers in Behavioral Neuroscience 4, 21 (2010). DOI: [10.3389/fnbeh.2010.00021](https://doi.org/10.3389/fnbeh.2010.00021)
3. J Neuroscience 30, 32 (2010), DOI: [10.1523/JNEUROSCI.1818-10.2010](https://doi.org/10.1523/JNEUROSCI.1818-10.2010)
4. J Neurophysiol 122, 4 (2019), DOI: [10.1152/jn.00038.2019](https://doi.org/10.1152/jn.00038.2019)

This datasheet was generated on January 16, 2026 at 03:31:52 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/tetrodotoxin/>