

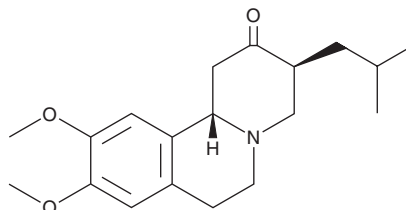
# PRODUCT INFORMATION



## Tetrabenazine

Item No. 20380

**CAS Registry No.:** 58-46-8  
**Formal Name:** *rel*-1,3R,4,6,7,11bR-hexahydro-9,10-dimethoxy-3-(2-methylpropyl)-2H-benzo[a]quinolizin-2-one  
**Synonyms:** NSC 169886, NSC 172187, Ro 1-9569, TBZ  
**MF:** C<sub>19</sub>H<sub>27</sub>NO<sub>3</sub>  
**FW:** 317.4  
**Purity:** ≥98% (mixture of enantiomers)  
**UV/Vis.:** λ<sub>max</sub>: 285 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Tetrabenazine (TBZ) is supplied as a crystalline solid. A stock solution may be made by dissolving the TBZ in the solvent of choice, which should be purged with an inert gas. TBZ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TBZ in these solvents is approximately 10, 25, and 30 mg/ml, respectively.

TBZ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TBZ should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TBZ has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Tetrabenazine is an inhibitor of vesicular monoamine transporter 2 (VMAT2) that is selective over VMAT1 (K<sub>s</sub> = 97 and >20,000 nM, respectively, in a serotonin uptake assay).<sup>1</sup> It dose-dependently reduces levels of the monoamines norepinephrine (Item No. 16673), dopamine (Item No. 21992), and serotonin (5-HT; Item No. 14332) in rat brain and has been used to induce depressive-like behavior in animal models.<sup>2,3</sup> Tetrabenazine (5 mg/kg) improves performance in balance beam and rotarod tests and prevents decreases in the number of striatal medium spiny neurons (MSNs) in a YAC128 transgenic mouse model of Huntington's disease.<sup>4</sup> Formulations containing tetrabenazine have been used in the treatment of chorea associated with Huntington's disease.

### References

1. Erickson, J.D., Schäfer, M.K.H., Bonner, T.I., *et al.* Distinct pharmacological properties and distribution in neurons and endocrine cells of two isoforms of the human vesicular monoamine transporter. *Proc. Natl. Acad. Sci. U.S.A.* **93**(10), 5166-5171 (1996).
2. Pettibone, D.J., Totaro, J.A., and Pflueger, A.B. Tetrabenazine-induced depletion of brain monoamines: Characterization and interaction with selected antidepressants. *Eur. J. Pharmacol.* **102**(3-4), 425-430 (1984).
3. Preskorn, S.H., Kent, T.A., Glotzbach, R.K., *et al.* Cerebromicrocirculatory defects in animal model of depression. *Psychopharmacology (Berl.)* **84**(2), 196-199 (1984).
4. Wang, H., Chen, X., Li, Y., *et al.* Tetrabenazine is neuroprotective in Huntington's disease mice. *Mol. Neurodegener.* **5**, 18 (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, 12/01/2022

#### 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](http://WWW.CAYMANCHEM.COM)