

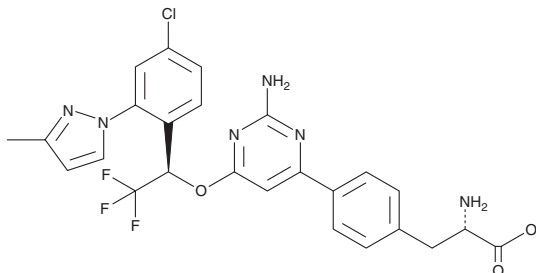
PRODUCT INFORMATION



Telotristat

Item No. 29424

CAS Registry No.: 1033805-28-5
Formal Name: 4-[2-amino-6-[(1R)-1-[4-chloro-2-(3-methyl-1H-pyrazol-1-yl)phenyl]-2,2,2-trifluoroethoxy]-4-pyrimidinyl]-L-phenylalanine
Synonym: LP778902
MF: C₂₅H₂₂ClF₃N₆O₃
FW: 546.9
Purity: ≥98%
UV/Vis.: λ_{max}: 246, 307 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

Telotristat is supplied as a crystalline solid. A stock solution may be made by dissolving the telotristat in the solvent of choice, which should be purged with an inert gas. Telotristat is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of telotristat in ethanol is approximately 0.5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Telotristat is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, telotristat should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Telotristat has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Telotristat is a tryptophan hydroxylase 1 (TPH1) inhibitor (IC₅₀ = 16 nM).¹ Dietary administration of telotristat to pregnant and lactating mice decreases serotonin (5-HT; Item No. 14332) levels in maternal serum and decreases the femoral bone volume to tissue volume ratio, trabecular number and thickness, and cortical tissue mineral density, as well as serum calcium levels, in pups.²

References

1. Goldberg, D.R., De Lombaert, S., Aiello, R., *et al.* Discovery of acyl guanidine tryptophan hydroxylase-1 inhibitors. *Bioorg. Med. Chem. Lett.* **26**(12), 2855-2860 (2016).
2. Weaver, S.R., Fricke, H.P., Xie, C., *et al.* Peripartum dietary supplementation of a small-molecule inhibitor of tryptophan hydroxylase 1 compromises infant, but not maternal, bone. *Am. J. Physiol. Endocrinol. Metab.* **215**(6), E1133-E1142 (2018).

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, 01/03/2023

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