

Telotristat

Catalog #: 27340

Lot #: 230317

Size: 5 mg

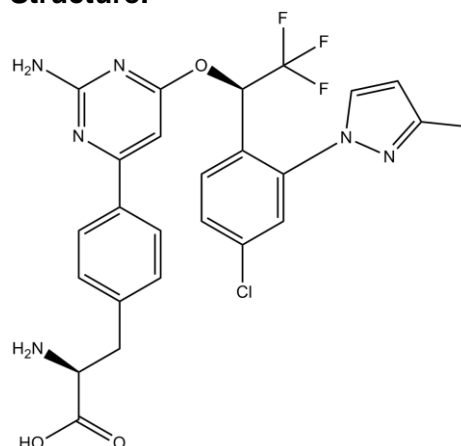
CAS Registry #: 1033805-28-5

Purity: ≥98%

Chemical Formula: C₂₅H₂₂ClF₃N₆O₃

Molecular Weight: 546.93 Da

Structure:



Description: Telotristat is the active metabolite of LX1606 (Telotristat etiprate), which is an orally bioavailable, small-molecule, tryptophan hydroxylase (TPH) inhibitor with potential antiserotonergic activity.

Synonym: LX1032

Appearance: White powder

Solubility: ≥27.35 mg/mL in DMSO with ultrasonic; insoluble in H₂O; insoluble in EtOH

Biological Activity: Telotristat has been reported to inhibit the TPH enzyme. This inhibition can reduce serotonin production in the periphery.

Storage/Stability: Store as supplied at or below -20°C for up to 2 years. Stable in anhydrous DMSO at -80°C for up to 6 months.

References:

1. Kulke, M., *et al.*, *Endocr. Relat. Cancer*, 2014; **21**: 705-714.
2. Margolis, K.G. *et al.*, *Gut*, 2014; **63(6)**:928-37.