

## Telotristat

**Catalog #:** 27340

**Lot #:** 160706

**Size:** 10 mg

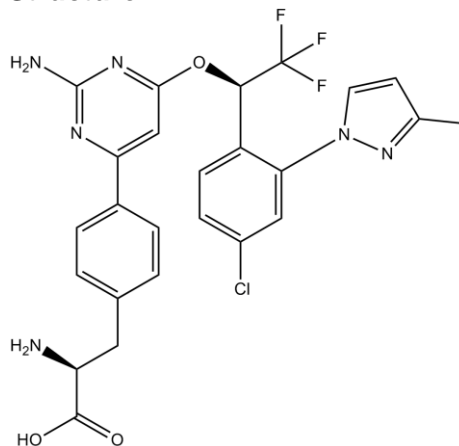
**CAS Registry #:** 1033805-28-5

**Structure:**

**Purity:** ≥98%

**Chemical Formula:** C<sub>25</sub>H<sub>22</sub>ClF<sub>3</sub>N<sub>6</sub>O<sub>3</sub>

**Molecular Weight:** 546.9



**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:** Soluble in DMSO (10 mM). Only sparingly soluble in water (<1.5 µg/ml).

**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.