

# PRODUCT INFORMATION

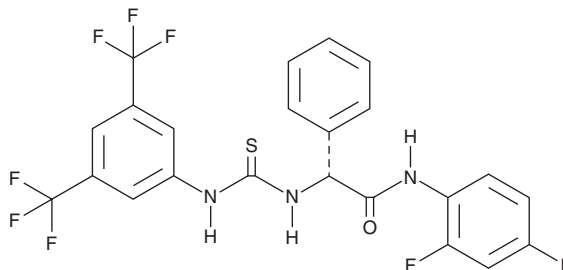


## BL 918

Item No. 34303

**CAS Registry No.:** 2101517-69-3  
**Formal Name:**  $\alpha$ R-[[[3,5-bis(trifluoromethyl)phenyl]amino]thioxomethyl]amino]-N-(2,4-difluorophenyl)-benzeneacetamide

**MF:** C<sub>23</sub>H<sub>15</sub>F<sub>8</sub>N<sub>3</sub>OS  
**FW:** 533.4  
**Purity:** ≥98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 251 nm  
**Supplied as:** A 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

BL 918 is supplied as a solid. A stock solution may be made by dissolving the BL 918 in the solvent of choice, which should be purged with an inert gas. BL 918 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BL 918 in these solvents is approximately 30 mg/ml.

### Description

BL 918 is an activator of unc-51-like autophagy activating kinase 1 (ULK1).<sup>1</sup> It activates ULK1 with an EC<sub>50</sub> value of 24.14 nM in an ADP-based kinase assay. BL 918 (5  $\mu$ M) increases the levels of ULK1 phosphorylated at serine 317 (Ser<sup>317</sup>) or Ser<sup>555</sup>, reduces the level of ULK1 phosphorylated at Ser<sup>757</sup>, and induces autophagy in SH-SY5Y cells. It reduces motor dysfunction and increases the number of dopaminergic neurons in the striatum in a mouse model of Parkinson's disease induced by MPTP when administered at doses of 40 and 80 mg/kg.

### Reference

1. Ouyang, L., Zhang, L., Zhang, S., *et al.* Small-molecule activator of UNC-51-like kinase 1 (ULK1) that induces cytoprotective autophagy for Parkinson's disease treatment. *J. Med. Chem.* **61(7)**, 2776-2792 (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, 12/08/2022

### 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