

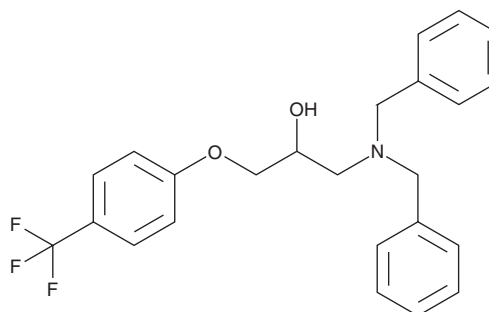
# PRODUCT INFORMATION



## BC-1618

Item No. 36136

**CAS Registry No.:** 2222094-18-8  
**Formal Name:** 1-[bis(phenylmethyl)amino]-3-[4-(trifluoromethyl)phenoxy]-2-propanol  
**MF:** C<sub>24</sub>H<sub>24</sub>F<sub>3</sub>NO<sub>2</sub>  
**FW:** 415.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 227 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

BC-1618 is supplied as a solid. A stock solution may be made by dissolving the BC-1618 in the solvent of choice, which should be purged with an inert gas. BC-1618 is soluble in DMSO.

### Description

BC-1618 is an inhibitor of F-box only protein 48 (FBXO48), an E3 ubiquitin ligase.<sup>1</sup> It is selective for FBXO48 over FBXO30 at 3 μM. BC-1618 (0.5-10 μM) increases the levels of phosphorylated AMP-activated protein kinase (pAMPK) and phosphorylated acetyl-CoA carboxylase (pACC), an AMPK downstream substrate, in HEK293T cells expressing human FBXO48. It induces mitochondrial fission in BEAS-2B cells and autophagy in HEK293A cells expressing GFP-tagged LC3. BC-1618 also inhibits LPS-induced activation of NF-κB in THP-1 cells in a reporter assay and cytokine release in isolated human peripheral blood mononuclear cells (PBMCs). It increases insulin sensitivity in a mouse model of high-fat diet-induced obesity when administered at a dose of 20 mg/kg.

### Reference

1. Liu, Y., Jurczak, M.J., Lear, T.B., *et al.* A Fbxo48 inhibitor prevents pAMPKα degradation and ameliorates insulin resistance. *Nat. Chem. Biol.* **17**(3), 298-306 (2021).

#### 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, 12/19/2022

#### 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