

# PRODUCT INFORMATION



## 84-B10

Item No. 43187

CAS Registry No.: 698346-43-9

Formal Name:  $\beta$ -[2-[[2-(4-methoxyphenoxy)-5-(trifluoromethyl)phenyl]amino]-2-oxoethyl]-benzenepropanoic acid

MF:  $C_{25}H_{22}F_3NO_5$

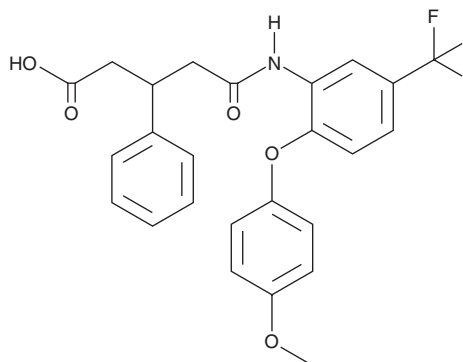
FW: 473.4

Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^\circ\text{C}$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

84-B10 is supplied as a solid. A stock solution may be made by dissolving the 84-B10 in the solvent of choice, which should be purged with an inert gas. 84-B10 is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

84-B10 is a derivative of 3-phenylglutaric acid.<sup>1</sup> It restores mitochondrial architecture and increases spare respiratory capacity in TKPTS renal tubular epithelial cells when used at concentrations ranging from 10 to 100  $\mu\text{M}$ . 84-B10 (10-100  $\mu\text{M}$ ) inhibits lipid peroxidation and ferroptosis induced by cisplatin (Item No. 13119) in the same cells. *In vivo*, 84-B10 (5, 10, and 15 mg/kg) reduces renal tubular injury and blood urea nitrogen (BUN) levels in a mouse model of cisplatin-induced acute kidney injury.

### Reference

1. Fan, J., Xu, X., Li, Y., *et al.* A novel 3-phenylglutaric acid derivative (84-B10) alleviates cisplatin-induced acute kidney injury by inhibiting mitochondrial oxidative stress-mediated ferroptosis. *Free. Radic. Biol. Med.* **194**, 84-98 (2023).

#### 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, 03/27/2025

#### 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