

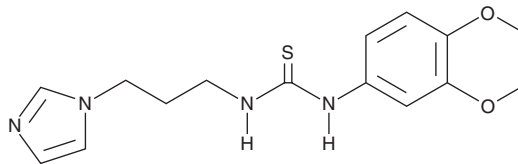
PRODUCT INFORMATION



PBD-150

Item No. 36112

CAS Registry No.: 790663-33-1
Formal Name: N-(3,4-dimethoxyphenyl)-N'-[3-(1H-imidazol-1-yl)propyl]-thiourea
MF: C₁₅H₂₀N₄O₂S
FW: 320.4
Purity: ≥95%
UV/Vis.: λ_{max}: 249 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

PBD-150 is supplied as a solid. A stock solution may be made by dissolving the PBD-150 in the solvent of choice, which should be purged with an inert gas. PBD-150 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PBD-150 in these solvents is approximately 1 mg/ml.

Description

PBD-150 is an inhibitor of glutaminyl cyclase ($K_i = 0.49 \mu\text{M}$ for the human enzyme).¹ It inhibits the formation of the pyroglutamic acid-containing amyloid- β (A β) peptides pGlu-A β (3-40) and pGlu-A β (3-42) in HEK293 cells expressing human glutaminyl cyclase and human amyloid precursor protein (APP) with an N-terminal deletion of the first two amino acids.¹

Reference

1. Buchholz, M., Heiser, U., Schilling, S., *et al.* The first potent inhibitors for human glutaminyl cyclase: Synthesis and structure-activity relationship. *J. Med. Chem.* **49**(2), 664-677 (2006).

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