

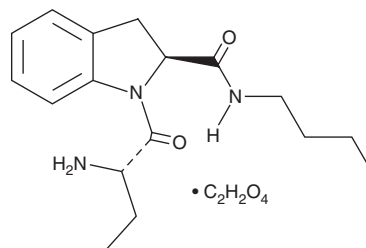
PRODUCT INFORMATION



Butabindide (oxalate)

Item No. 21610

CAS Registry No.: 185213-03-0
Formal Name: [S-(R*,R*)]-1-(2-amino-1-oxobutyl)-N-butyl-2,3-dihydro-1H-indole-2-carboxamide, monoethanedioate
MF: C₁₇H₂₅N₃O₂ • C₂H₂O₄
FW: 393.4
Purity: ≥98%
UV/Vis.: λ_{max}: 250, 280, 288 nm
Supplied as: A crystalline 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

Butabindide (oxalate) is supplied as a crystalline solid. A stock solution may be made by dissolving the butabindide (oxalate) in the solvent of choice, which should be purged with an inert gas. Butabindide (oxalate) is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 1.6 mg/ml.

Butabindide (oxalate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, butabindide (oxalate) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Butabindide (oxalate) has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Butabindide is a potent inhibitor of cholecystokinin-inactivating peptidase/tripeptidyl peptidase 2 (CCK-inactivating peptidase/TPP-2; K_i = 7 nM).¹ It is selective for CCK-inactivating peptidase/TPP-2 over a panel of serine proteases (K_is = >1 μM) as well as CCK receptors (K_is = >0.1 mM). Butabindide (0.1-100 μM) increases levels of CCK octapeptide (Item Nos. 24404 | 23371) in depolarized rat cerebral cortex slices. *In vivo*, butabindide inhibits CCK-inactivating peptidase/TPP-2 in mouse liver and brain (ID₅₀s = 1.1 and 6.8 mg/kg, respectively). It also enhances CCK octapeptide-induced delay in gastric emptying and reduces food intake in mice.

Reference

1. Rose, C., Vargas, F., Facchinetti, P., *et al.* Characterization and inhibition of a cholecystokinin-inactivating serine peptidase. *Nature* **380(6573)**, 403-409 (1996).

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, 10/06/2022

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