

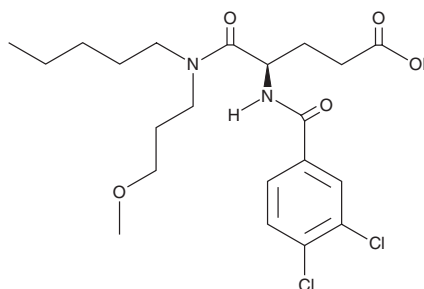
PRODUCT INFORMATION



Dexloxiglumide

Item No. 34759

CAS Registry No.: 119817-90-2
Formal Name: (4R)-4-[(3,4-dichlorobenzoyl)amino]-5-[(3-methoxypropyl)pentylamino]-5-oxo-pentanoic acid
MF: C₂₁H₃₀Cl₂N₂O₅
FW: 461.4
Purity: ≥98%
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

Dexloxiglumide is supplied as a solid. A stock solution may be made by dissolving the dexloxiglumide in the solvent of choice, which should be purged with an inert gas. Dexloxiglumide is soluble in the organic solvent DMSO (sonicated) at a concentration of approximately 50 mg/ml.

Description

Dexloxiglumide is a cholecystokinin (CCK) receptor antagonist and is the single (R) isomer of loxiglumide (Item No. 25534).^{1,2} It selectively binds to the CCK₁ receptor over CCK₂ in CHO cells expressing the rat receptors (K_s = 0.0234 and 5.88 μM, respectively).¹ Dexloxiglumide (20 μmol/kg) inhibits guinea pig gallbladder contractions induced by the CCK agonist CCK octapeptide (sulfated) (CCK-8S; Item No. 23371). It reduces CCK-8-induced gastric emptying in rats (ID₅₀ = 1.14 mg/kg).²

References

1. Morton, M.F., Barrett, T.D., Yan, W., *et al.* 3-[5-(3,4-Dichloro-phenyl)-1-(4-methoxy-phenyl)-1H-pyrazol-3-yl]-2-m-tolyl-propionate (JNJ-17156516), a novel, potent, and selective cholecystokinin 1 receptor antagonist: In vitro and in vivo pharmacological comparison with dexloxiglumide. *J. Pharmacol. Exp. Ther.* **323**(2), 562-569 (2007).
2. Scarpignato, C., Kisfalvi, I., D'Amato, M., *et al.* Effect of dexloxiglumide and spiroglumide, two new CCK-receptor antagonists, on gastric emptying and secretion in the rat: Evaluation of their receptor selectivity *in vivo*. *Aliment. Pharmacol. Ther.* **10**(3), 411-419 (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, 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