

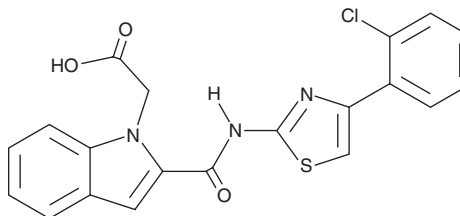
PRODUCT INFORMATION



SR 27897

Item No. 28511

CAS Registry No.: 136381-85-6
Formal Name: 2-[[[4-(2-chlorophenyl)-2-thiazolyl]amino]carbonyl]-1H-indole-1-acetic acid
Synonym: Lintitript
MF: C₂₀H₁₄ClN₃O₃S
FW: 411.9
Purity: ≥98%
UV/Vis.: λ_{max}: 321 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

SR 27897 is supplied as a solid. A stock solution may be made by dissolving the SR 27897 in the solvent of choice, which should be purged with an inert gas. SR 27897 is soluble in DMSO.

Description

SR 27897 is a nonpeptide cholecystokinin (CCK) receptor antagonist.¹ It selectively binds to CCK₁ receptors in rat pancreatic membranes (IC₅₀ = 0.58 nM) over CCK₂ receptors in guinea pig cortical membranes and gastrin receptors in guinea pig gastric gland suspensions (IC₅₀s = 489 and 2,883 nM, respectively). SR 27897 inhibits amylase secretion induced by CCK in isolated rat pancreatic acini (pA₂ = 7.50) and CCK-induced guinea pig gallbladder contractions *ex vivo* (pA₂ = 9.57). It completely reverses CCK-induced amylase secretion in rats when administered at a dose of 1 mg/kg. SR 27897 also inhibits CCK-induced gastric and gallbladder emptying in mice (ED₅₀s = 3 and 72 µg/kg, respectively). SR 27897 inhibits gallbladder emptying in a mouse model of egg yolk-stimulated endogenous CCK release (ED₅₀ = 27 µg/kg).

Reference

1. Gully, D., Fréhel, D., Marcy, C., et al. Peripheral biological activity of SR 27897: A new potent non-peptide antagonist of CCKA receptors. *Eur. J. Pharmacol.* **232**(1), 13-19 (1993).

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, 11/22/2022

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