

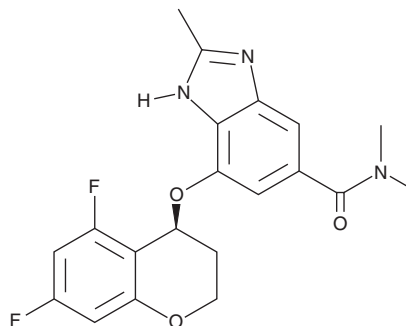
PRODUCT INFORMATION



Tegoprazan

Item No. 33243

CAS Registry No.: 942195-55-3
Formal Name: 7-[[[(4S)-5,7-difluoro-3,4-dihydro-2H-1-benzopyran-4-yl]oxy]-N,N,2-trimethyl-1H-benzimidazole-5-carboxamide
Synonym: CJ-12420
MF: $C_{20}H_{19}F_2N_3O_3$
FW: 387.4
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 220 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

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

Description

Tegoprazan is a proton pump inhibitor that inhibits H^+/K^+ -ATPase activity in porcine gastric membrane vesicles ($\text{IC}_{50} = 0.53 \mu\text{M}$).¹ It completely inhibits basal gastric acid secretion in pylorus-ligated rats when administered at a dose of 10 mg/kg and reduces esophageal mucosal tissue levels of myeloperoxidase (MPO), a marker of tissue damage, in a rat model of reflux esophagitis ($\text{ED}_{50} = 5.9 \text{ mg/kg}$). Tegoprazan reduces ulcer area in rat models of gastric ulcers induced by naproxen, water-immersion restraint stress (WIRS), or ethanol (ED_{50} s = 0.1, 0.1, and 1.4 mg/kg, respectively).

Reference

1. Kim, D.K., Lee, K.-H., Kim, S.-J., *et al.* Effects of tegoprazan, a novel potassium-competitive acid blocker, on rat models of gastric acid-related disease. *J. Pharmacol. Exp. Ther.* **369**(3), 318-327 (2019).

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, 11/14/2022

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