

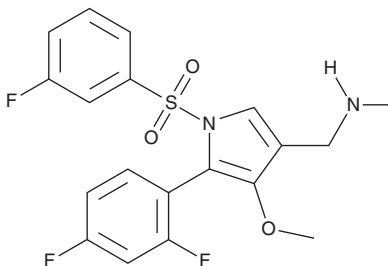
PRODUCT INFORMATION



Abeprazan

Item No. 45283

CAS Registry No.: 1902954-60-2
Formal Name: 5-(2,4-difluorophenyl)-1-[(3-fluorophenyl)sulfonyl]-4-methoxy-N-methyl-1H-pyrrole-3-methanamine
Synonyms: DWP14012, Fexuprazan
MF: C₁₉H₁₇F₃N₂O₃S
FW: 410.4
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

Abeprazan is supplied as a solid. A stock solution may be made by dissolving the abeprazan in the solvent of choice, which should be purged with an inert gas. Abeprazan is slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water and DMSO.

Description

Abeprazan is a potassium-competitive acid blocker (P-CAB).^{1,2} Abeprazan (20 µg/ml) increases *trans*-epithelial electrical resistance (TEER) and levels of the tight junction proteins occludin and mucin-1 (Muc-1) in a Caco-2 colon cell model of ulcerative colitis (UC) induced by dextran sulfate sodium (DSS; Item No. 23250).¹ *In vivo*, abeprazan (20 mg/kg) reduces ileal levels of TNF-α, IL-6, and IL-1β in a mouse model of small intestinal injury induced by indomethacin (IDM; Item No. 70270).² Formulations containing abeprazan have been used in the treatment of erosive esophagitis (EE) and gastroesophageal reflux disease (GERD).

References

1. Jeon, J.-H., Lee, J., Baek, Y.-J., *et al.* Restoration of intestinal barrier function by fexuprazan, a potassium-competitive acid blocker, in Caco-2 cells and its higher gastrointestinal distribution in rats. *Biomed. Pharmacother.* **188**, 118185 (2025).
2. Lee, Y., Kim, Y., Park, S., *et al.* Fexuprazan mitigates NSAID-induced small intestinal injury by restoring intestinal barrier integrity in mice. *Biomed. Pharmacother.* **190**, 118386 (2025).

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, 02/25/2026

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