

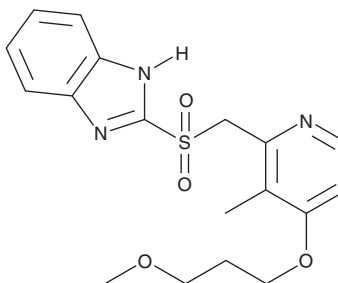
# PRODUCT INFORMATION



## Rabeprazole sulfone

Item No. 35076

CAS Registry No.: 117976-47-3  
Formal Name: 2-[[[4-(3-methoxypropoxy)-3-methyl-2-pyridinyl]methyl]sulfonyl]-1H-benzimidazole  
MF:  $C_{18}H_{21}N_3O_4S$   
FW: 375.4  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Rabeprazole sulfone is supplied as a solid. A stock solution may be made by dissolving the rabeprazole sulfone in the solvent of choice, which should be purged with an inert gas. Rabeprazole sulfone is slightly soluble in methanol and acetone.

### Description

Rabeprazole sulfone is a metabolite of the proton pump inhibitor rabeprazole (Item No. 14939).<sup>1</sup> It is formed from rabeprazole by the cytochrome P450 (CYP) isoform CYP3A. Rabeprazole sulfone has also been found as an impurity in bulk preparations of rabeprazole.<sup>2</sup>

### References

1. VandenBranden, M., Ring, B.J., Binkley, S.N., *et al.* Interaction of human liver cytochromes P450 *in vitro* with LY307640, a gastric proton pump inhibitor. *Pharmacogenetics* **6**(1), 81-91 (1996).
2. Reddy, G.M., Bhaskar, B.V., Reddy, P.P., *et al.* Identification and characterization of potential impurities of rabeprazole sodium. *J. Pharm. Biomed. Anal.* **43**(4), 1262-1269 (2007).

#### 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, 12/09/2022

#### 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