

PRODUCT INFORMATION



Sucralfate (technical grade)

Item No. 27882

CAS Registry No.: 54182-58-0

Formal Name: hexadeca- μ -hydroxytetracosahydroxy[μ_8 -[[1,3,4,6-tetra-O-sulfo- β -D-fructofuranosyl α -D-glucopyranoside tetrakis(sulfato- κ O)](8-)]hexadeca-aluminum

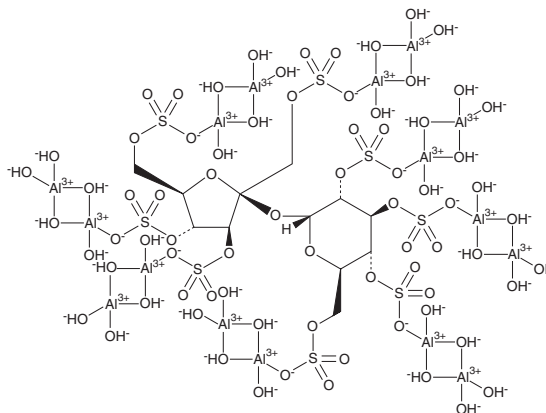
MF: $C_{12}H_{54}Al_{16}O_{75}S_8$

FW: 2,086.7

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

Sucralfate (technical grade) is supplied as a solid. A stock solution may be made by dissolving the sucralfate (technical grade) in the solvent of choice, which should be purged with an inert gas. Sucralfate (technical grade) is slightly soluble in the aqueous acids.

Description

Sucralfate is a basic aluminum sucrose sulfate complex that has gastroprotective activity.¹ It inhibits rat pepsin in a concentration-dependent manner and pepsin activity in isolated human gastric juice. It also inhibits ulcer formation induced by pyloric ligation, indomethacin (Item No. 70270), or cysteamine (Item No. 22193) in rats. Sucralfate (5,600 mg/animal) is protective against neutral ethanol and acidified taurocholic acid-induced damage in a rat model of hydrochloric acid-induced gastric mucosal damage, increasing the pH and reducing the disappearance of hydrogen ions.² Formulations containing sucralfate have been used as antacids in the treatment of duodenal ulcer.

References

1. Borella, L.E., Seethaler, K., and Lippmann, W. Sucralfate: Antipeptic, antiulcer activities and antagonism of gastric emptying. *Arzneimittelforschung* **29(5)**, 793-798 (1979).
2. Harrington, S.J., Schlegel, J.F., and Code, C.F. The protective effect of sucralfate on the gastric mucosa of rats. *J. Clin. Gastroenterol.* **3(Suppl 2)**, 129-134 (1981).

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, 01/17/2023

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