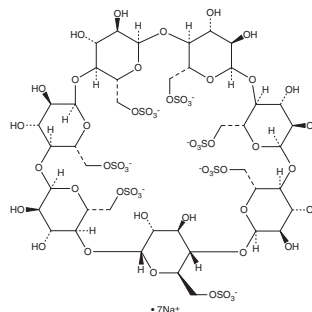


PRODUCT INFORMATION

6-O-Sulfo-β-cyclodextrin (sodium salt)

Item No. 30168

CAS Registry No.: 197587-31-8
Formal Name: 6^A,6^B,6^C,6^D,6^E,6^F,6^G-heptakis(hydrogen sulfate), β-cyclodextrin, heptasodium salt
Synonyms: Heptakis(6-O-sulfo)-β-cyclodextrin, HS-β-CD
MF: C₄₂H₆₃O₅₆S₇ • 7Na
FW: 1,849.3
Purity: ≥95%
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

6-O-Sulfo-β-cyclodextrin (sodium salt) is supplied as a solid. Aqueous solutions of 6-O-sulfo-β-cyclodextrin (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 6-O-sulfo-β-cyclodextrin (sodium salt) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

6-O-Sulfo-β-cyclodextrin is a sulfated cyclodextrin.¹ It has been used for the chiral separation of nadolol racemates by capillary electrophoresis.

Reference

1. Wang, F., Dowling, T., Bicker, G., *et al.* Electrophoretic chiral separation of pharmaceutical compounds with multiple stereogenic centers in charged cyclodextrin media. *J. Sep. Sci.* **24**(5), 378-384 (2001).

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, 12/13/2022

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