

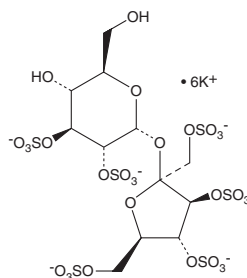
# PRODUCT INFORMATION



## Sucrose hexasulfate (potassium salt)

Item No. 18239

**CAS Registry No.:** 386229-70-5  
**Formal Name:** 1,3,4,6-tetra-O-sulfo-β-D-fructofuranosyl α-D-glucopyranoside 2,3-bis(hydrogen sulfate), hexapotassium salt  
**MF:** C<sub>12</sub>H<sub>16</sub>O<sub>29</sub>S<sub>6</sub> • 6K  
**FW:** 1,051.2  
**Purity:** ≥65%  
**Supplied as:** A crystalline 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

Sucrose hexasulfate (potassium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the sucrose hexasulfate (potassium salt) in water. The solubility of sucrose hexasulfate (potassium salt) in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Sucralfate is a gastrointestinal protectant that includes sucrose octasulfate (Item No. 16382). Sucralfate protects gastric epithelial cells against acid- and pepsin-induced damage.<sup>1</sup> Sucrose hexasulfate is a polysulfated disaccharide that is used as a reference standard for sucralfate.<sup>2</sup>

### References

1. Furukawa, O., Matsui, H., and Suzuki, N. Effects of sucralfate and its components on acid- and pepsin-induced damage to rat gastric epithelial cells. *Jpn. J. Pharmacol.* **75(1)**, 21-25 (1997).
2. Silvey, G.L. Two-dimensional and deuterium-induced, differential-isotope-shift nuclear magnetic resonance characterization of potassium sucrose octasulfate and sucralfate. *J. Pharm. Sci.* **81(5)**, 471-474 (1992).

#### 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, 11/10/2022

#### 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