

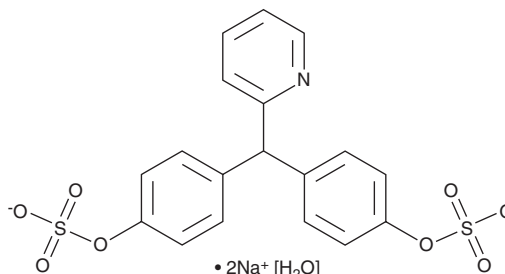
# PRODUCT INFORMATION



## Picosulfate (sodium salt hydrate)

Item No. 34566

**CAS Registry No.:** 1307301-38-7  
**Formal Name:** 4,4'-(2-pyridinylmethylene)bis-phenol, 1,1'-bis(hydrogen sulfate), disodium salt, monohydrate  
**MF:** C<sub>18</sub>H<sub>13</sub>NO<sub>8</sub>S<sub>2</sub> • 2Na [H<sub>2</sub>O]  
**FW:** 499.4  
**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

Picosulfate (sodium salt hydrate) is supplied as a solid. A stock solution may be made by dissolving the picosulfate (sodium salt hydrate) in the solvent of choice, which should be purged with an inert gas. Picosulfate (sodium salt hydrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of picosulfate (sodium salt hydrate) in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of picosulfate (sodium salt hydrate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of picosulfate (sodium salt hydrate) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Picosulfate is a stimulant laxative and prodrug form of desacetyl bisacodyl (DAB; Item No. 20928).<sup>1</sup> Unlike bisacodyl (Item No. 28826), picosulfate is hydrolyzed to DAB by colonic microbiota. Picosulfate (5 mg/kg) increases excreted fecal weight in a rat model of loperamide-induced constipation.<sup>2</sup> Formulations containing picosulfate have been used for colonoscopy preparation.

### References

1. Twycross, R., Sykes, N., Mihalyo, M., *et al.* Stimulant laxatives and opioid-induced constipation. *J. Pain Symptom Manage.* **43**(2), 306-313 (2012).
2. Méité, S., Bahi, C., Yéo, D., *et al.* Laxative activities of *Mareya micrantha* (Benth.) Müll. Arg. (Euphorbiaceae) leaf aqueous extract in rats. *BMC Complement. Altern. Med.* **10**, 7 (2010).

#### 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, 11/04/2022

#### 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