

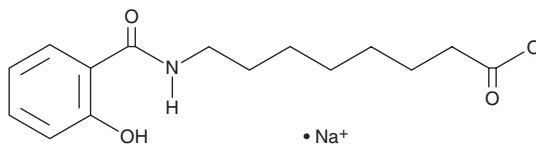
# PRODUCT INFORMATION



## Salcaprozate (sodium salt)

Item No. 36505

**CAS Registry No.:** 203787-91-1  
**Formal Name:** 8-[(2-hydroxybenzoyl)amino]-octanoic acid, monosodium salt  
**Synonym:** SNAC  
**MF:** C<sub>15</sub>H<sub>20</sub>NO<sub>4</sub> • Na  
**FW:** 301.3  
**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

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

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 salcaprozate (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of salcaprozate (sodium salt) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Salcaprozate is an intestinal permeation enhancer.<sup>1</sup> It increases jejunal accumulation of the bioactive antioxidant peptide PW5 *in vitro* in a non-everted rat gut sac model. *In vivo*, salcaprozate increases plasma concentrations of orally administered human parathyroid hormone (1-34) in pigs.<sup>2</sup>

### References

- Xu, Y., Amakye, W.K., Xiao, G., *et al.* Intestinal absorptivity-increasing effects of sodium N-[8-(2-hydroxybenzoyl)amino]-caprylate on food-derived bioactive peptide *Food Chem.* **401**, 134059 (2023).
- Burshtein, G., Itin, C., Tang, J.C.Y., *et al.* The combined effect of permeation enhancement and proteolysis inhibition on the systemic exposure of orally administered peptides: Salcaprozate sodium, soybean trypsin inhibitor, and teriparatide study in pigs. *Int. J. Pharm. X.* **3**, 100097 (2021).

#### 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, 05/02/2023

#### 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