

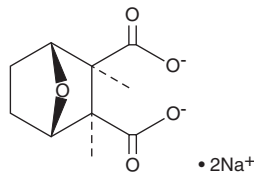
# PRODUCT INFORMATION



## Cantharidic Acid (sodium salt)

Item No. 19073

**CAS Registry No.:** 1465-77-6  
**Formal Name:** (1R,2S,3R,4S)-rel-2,3-dimethyl-7-oxabicyclo[2.2.1]heptane-2,3-dicarboxylic acid, disodium salt  
**MF:** C<sub>10</sub>H<sub>12</sub>O<sub>5</sub> • 2Na  
**FW:** 258.2  
**Purity:** ≥95%  
**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

Cantharidic acid (sodium salt) is supplied as a crystalline solid. Aqueous solutions of cantharidic acid (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of cantharidic acid (sodium salt) in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Cantharidic acid is a hydrolysis product of cantharidin (Item No. 14589), an inhibitor of protein phosphatases first isolated from Chinese blister beetles.<sup>1</sup> Cantharidic acid inhibits the protein phosphatases PP1 and PP2A with IC<sub>50</sub> values of 0.6 and 0.05 μM, respectively, but does not affect PP2B or PP2C.<sup>1</sup> Cantharidic acid is frequently used to study cellular processes that are regulated by reversible protein phosphorylation.

### Reference

1. McCluskey, A., Sim, A.T.R., and Sakoff, J.A. Serine-threonine protein phosphatase inhibitors: Development of potential therapeutic strategies. *J. Med. Chem.* **45(6)**, 1151-1175 (2002).

#### 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, 03/15/2023

#### 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