

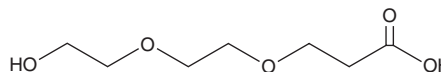
# PRODUCT INFORMATION



## Hydroxy-PEG<sub>2</sub> Acid

Item No. 30903

CAS Registry No.: 1334286-77-9  
Formal Name: 3-[2-(2-hydroxyethoxy)ethoxy]-  
propanoic acid  
MF: C<sub>7</sub>H<sub>14</sub>O<sub>5</sub>  
FW: 178.2  
Purity: ≥95%  
Supplied as: A liquid  
Storage: -20°C  
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Hydroxy-PEG<sub>2</sub> acid is supplied as a liquid. A stock solution may be made by dissolving the hydroxy-PEG<sub>2</sub> acid in the solvent of choice, which should be purged with an inert gas. Hydroxy-PEG<sub>2</sub> acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Hydroxy-PEG<sub>2</sub> acid in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Hydroxy-PEG<sub>2</sub> acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, hydroxy-PEG<sub>2</sub> acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Hydroxy-PEG<sub>2</sub> acid has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Hydroxy-PEG<sub>2</sub> acid is a PEGylated building block.<sup>1,2</sup> It has been used in the synthesis of lysophosphatidic acid (LPA) receptor antagonists and ionic complexes.

### References

1. Sharma, S., Van Der Ploeg, L., and Henderson, B. Pharmaceutical compositions. *Rythem Metabolic, Inc.* **WO2014/144842 A2** (2014).
2. Bolli, M., Brotschi, C., Lescop, C., *et al.* Alkoxy-substituted pyridinyl derivatives as LPA1 receptor antagonists and their use in the treatment of fibrosis. *Idorsia Pharmaceuticals Ltd.* **WO2019/234115 A1** (2019).

#### 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, 04/18/2022

#### 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