

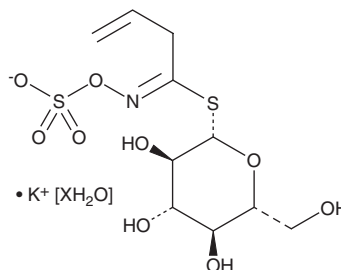
# PRODUCT INFORMATION



## Sinigrin (hydrate)

Item No. 26028

**Formal Name:** 1-[N-(sulfooxy)-3-butenimide]1-thio-β-D-glucopyranose, monopotassium salt, hydrate  
**MF:** C<sub>10</sub>H<sub>16</sub>NO<sub>9</sub>S<sub>2</sub> • K [XH<sub>2</sub>O]  
**FW:** 397.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 227 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Brassica juncea*



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

### Laboratory Procedures

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

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

### Description

Sinigrin is a glucosinolate that has been found in *Brassica* and has diverse biological activities.<sup>1-3</sup> It inhibits NOD-like receptor protein 3 (NLRP3) inflammasome activation in ATP-stimulated, LPS-primed RAW 264.7 cells in a concentration-dependent manner.<sup>1</sup> Sinigrin (1-100 μg/ml) reduces LPS-induced production of nitric oxide (NO) and prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010), as well as NF-κB activation, in RAW 264.7 cells. It induces cell cycle arrest at the G<sub>0</sub>/G<sub>1</sub> phase in HepG2 hepatocellular carcinoma cells when used at concentrations of 0.1 and 0.5 mM.<sup>2</sup> Sinigrin (10 and 20 mg/kg) reduces increases in renal glomerular basement membrane thickness, as well as increases in systolic and diastolic blood pressure, in a rat model of hypertension induced by angiotensin II.<sup>3</sup>

### References

1. Lee, H.-W., Lee, C.G., Rhee, D.-K., *et al.* Sinigrin inhibits production of inflammatory mediators by suppressing NF-κB/MAPK pathways or NLRP3 inflammasome activation in macrophages. *Int. Immunopharmacol.* **45**, 163-173 (2017).
2. Jie, M., Cheung, W.M., Yu, V., *et al.* Anti-proliferative activities of sinigrin on carcinogen-induced hepatotoxicity in rats. *PLoS One* **9(10)**, e110145 (2014).
3. Cong, C., Yuan, X., Hu, Y., *et al.* Sinigrin attenuates angiotensin II-induced kidney injury by inactivating nuclear factor-κB and extracellular signal-regulated kinase signaling *in vivo* and *in vitro*. *Int. J. Mol. Med.* **48(2)**, 161 (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, 01/03/2023

#### 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