

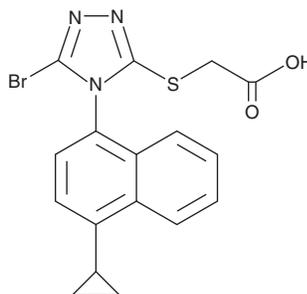
# PRODUCT INFORMATION



## Lesinurad

Item No. 23688

**CAS Registry No.:** 878672-00-5  
**Formal Name:** 2-[[5-bromo-4-(4-cyclopropyl-1-naphthalenyl)-4H-1,2,4-triazol-3-yl]thio]-acetic acid  
**MF:** C<sub>17</sub>H<sub>14</sub>BrN<sub>3</sub>O<sub>2</sub>S  
**FW:** 404.3  
**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

Lesinurad is supplied as a solid. A stock solution may be made by dissolving the lesinurad in the solvent of choice, which should be purged with an inert gas. Lesinurad is slightly soluble in organic solvents such as DMSO and methanol.

### Description

Lesinurad is an inhibitor of urate anion transporter 1 (URAT1; IC<sub>50</sub>s = 3.5 and 81 μM for human and rat transporters, respectively), a transporter that prevents renal urate resorption.<sup>1</sup> It decreases uptake of urate by MDCK cells transfected with human URAT1.<sup>2</sup> Lesinurad (80 mg/kg) also decreases serum levels of uric acid and urea nitrogen in a rat model of gout. Formulations containing lesinurad have been used in combination with other compounds in the treatment of gout.

### References

1. Tan, P.K., Ostertag, T.M., and Miner, J.N. Mechanism of high affinity inhibition of the human urate transporter URAT1. *Sci. Rep.* **6:34995**, (2016).
2. Wu, T., Chen, J., Dong, S., *et al.* Identification and characterization of a potent and selective inhibitor of human urate transporter 1. *Pharmacol. Rep.* (2017).

#### 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, 11/08/2022

#### 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