

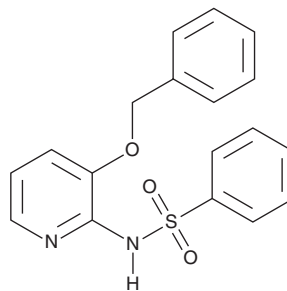
# PRODUCT INFORMATION



## TGN 073

Item No. 42086

**CAS Registry No.:** 877459-36-4  
**Formal Name:** N-[3-(phenylmethoxy)-2-pyridinyl]-benzenesulfonamide  
**MF:** C<sub>18</sub>H<sub>16</sub>N<sub>2</sub>O<sub>3</sub>S  
**FW:** 340.4  
**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

TGN 073 is supplied as a solid. A stock solution may be made by dissolving the TGN 073 in the solvent of choice, which should be purged with an inert gas. TGN 073 is soluble (≥10 mg/ml) in DMSO.

### Description

TGN 073 is a facilitator of aquaporin-4 (AQP4), an astroglial water channel.<sup>1,2</sup> It increases glymphatic transport and parenchymal penetration of a gadolinium-based tracer in rats when administered intraperitoneally at a dose of 200 mg/kg.<sup>1</sup> TGN 073 (200 mg/kg, i.p.) increases AQP4 polarization and reduces brain retention of exogenous amyloid-β (1-40) (Aβ40) in mice fed a high-fat diet.<sup>2</sup>

### References

1. Alghanimy, A., Martin, C., Gallagher, L., *et al.* The effect of a novel AQP4 facilitator, TGN-073, on glymphatic transport captured by diffusion MRI and DCE-MRI. *PLoS One* **18(3)**, e0282955 (2023).
2. Peng, W., Yuan, Y., Lei, J., *et al.* Long-term high-fat diet impairs AQP4-mediated glymphatic clearance of amyloid beta. *Mol. Neurobiol.* (2024).

#### 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, 11/27/2024

#### 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](http://WWW.CAYMANCHEM.COM)