

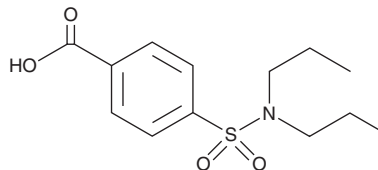
# PRODUCT INFORMATION



## Probenecid

Item No. 14981

CAS Registry No.: 57-66-9  
Formal Name: 4-[(dipropylamino)sulfonyl]-benzoic acid  
Synonym: NSC 18786  
MF:  $C_{13}H_{19}NO_4S$   
FW: 285.4  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{max}$ : 224, 249 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Probenecid is supplied as a crystalline solid. A stock solution may be made by dissolving the probenecid in the solvent of choice, which should be purged with an inert gas. Probenecid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of probenecid in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Probenecid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, probenecid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Probenecid has a solubility of approximately 0.15 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Probenecid is a benzoic acid derivative that inhibits organic anion transporters (OATs) and the ATP-binding cassette transporter (ABCC1) multidrug resistance protein 1 (MRP1) and activates the transient receptor potential (TRP) channel TRPV2.<sup>1-6</sup> It inhibits OAT1, OAT3, and OAT6 ( $K_s = 6.3, 9.0,$  and  $8.4 \mu M$ , respectively), as well as OAT2 ( $IC_{50} = 0.67 \mu M$ ), and is selective for OATs over organic cation transporter 1 (OCT1) and OCT2 ( $IC_{50}s = 1,600$  and  $1,700 \mu M$ , respectively).<sup>1-4</sup> Probenecid (7 mM) decreases Fas antibody-induced glutathione (GSH) efflux and staurosporine-induced efflux of the MRP1 substrate calcein in Jurkat cells and decreases apoptotic progression.<sup>6</sup> It is an agonist of TRPV2 ( $EC_{50} = 31.9 \mu M$ ).<sup>5</sup> It induces nociceptive behavior in mouse models of carrageenan- or complete Freund's adjuvant-induced inflammatory pain when administered at a dose of 20 mM in drinking water in combination with doses of carrageenan or CFA that do not provoke pain responses alone.<sup>5</sup> Formulations containing probenecid have been used in the treatment of gouty arthritis.

### References

1. Kaler, G., Truong, D.M., Khandelwal, A., et al. *J. Biol. Chem.* **282**(33), 23841-23853 (2007).
2. Takeda, M., Narikawa, S., Hosoyamada, M., et al. *Eur. J. Pharmacol.* **419**(2-3), 113-120 (2001).
3. Khamdang, S., Takeda, M., Shimoda, M., et al. *J. Pharmacol. Sci.* **94**(2), 197-202 (2004).
4. Arndt, P., Volk, C., Gorboulev, V., et al. *Am. J. Physiol. Renal Physiol.* **281**(3), F454-F468 (2001).
5. Bang, S., Kim, K.Y., Yoo, S., et al. *Neurosci. Lett.* **425**(2), 120-125 (2007).
6. Hammonnd, C.L., Marchan, R., Krance, S.M., et al. *J. Biol. Chem.* **282**(19), 14337-14347 (2007).

#### 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, 11/18/2022

#### 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