

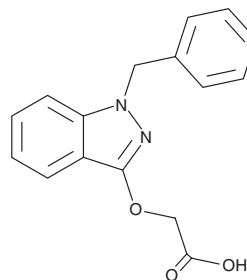
# PRODUCT INFORMATION



## Bendazac

Item No. 35127

**CAS Registry No.:** 20187-55-7  
**Formal Name:** 2-[[1-(phenylmethyl)-1H-indazol-3-yl]oxy]-acetic acid  
**Synonym:** AF 983  
**MF:** C<sub>16</sub>H<sub>14</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 282.3  
**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

Bendazac is supplied as a solid. A stock solution may be made by dissolving the bendazac in the solvent of choice, which should be purged with an inert gas. Bendazac is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of bendazac in these solvents is approximately 1, 15, and 20 mg/ml, respectively.

Bendazac is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, bendazac should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Bendazac has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Bendazac is a non-steroidal anti-inflammatory drug (NSAID).<sup>1,2</sup> It prevents denaturation of lens and serum proteins *in vitro*.<sup>1</sup> Bendazac (10-40 mM) prevents cyanate-induced opacification of the eye lens in a rat model of cataract.<sup>3</sup>

### References

1. Harding, J.J. Pharmacological treatment strategies in age-related cataracts. *Drugs Aging* **2(4)**, 287-300 (1992).
2. Guglielmotti, A., Capezone De Joannon, A., Cazzolla, N., *et al.* Radical scavenger activity of bendazac, an anticataract non-steroidal anti-inflammatory agent. *Pharmacol. Res.* **32(6)**, 369-373 (1995).
3. Lewis, B.S. and Harding, J.J. Evaluation of bendazac, a putative anti-cataract drug, in preventing the effects of cyanate in models of cataract. *T. Aging Res.* **6**, 127-130 (1986).

#### 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, 12/09/2022

#### 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