

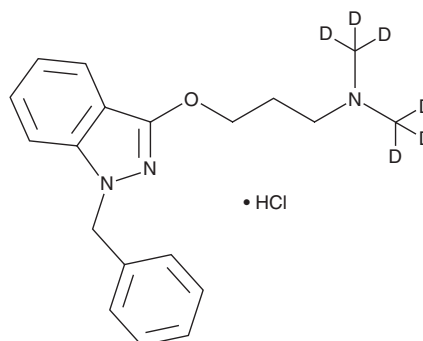
PRODUCT INFORMATION



Benzydamine-d₆ (hydrochloride)

Item No. 40112

CAS Registry No.: 1246817-20-8
Formal Name: N,N-dimethyl-3-[[1-(phenylmethyl-d₆)-1H-indazol-3-yl]oxy]-1-propanamine, monohydrochloride
MF: C₁₉H₁₇D₆N₃O • HCl
FW: 351.9
Chemical Purity: ≥98% (Benzydamine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₆); ≤1% d₀
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

Benzydamine-d₆ (hydrochloride) is intended for use as an internal standard for the quantification of benzydamine by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Benzydamine-d₆ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the benzydamine-d₆ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Benzydamine-d₆ (hydrochloride) is slightly soluble in chloroform and methanol.

Description

Benzydamine is a non-steroidal anti-inflammatory drug (NSAID).¹ It inhibits LPS-induced TNF-α and IL-1β production in isolated human peripheral blood mononuclear cells (PBMCs) when used at concentrations of 12, 25, and 50 μM. *In vivo*, benzydamine (20 mg/kg) reduces serum and lung levels of TNF-α and IL-1β, as well as mortality, in a mouse model of LPS-induced endotoxemia.² Benzydamine (100 mg/kg) reduces acetic acid-induced writhing in rats.³ Formulations containing benzydamine have been used as analgesics.

References

1. Sironi, M., Massimiliano, L., Transidico, P., *et al.* Differential effect of benzydamine on pro- versus anti-inflammatory cytokine production: Lack of inhibition of interleukin-10 and interleukin-1 receptor antagonist. *Int. J. Clin. Lab. Res.* **30**(1), 17-19 (2000).
2. Guglielmotti, A., Aquilini, L., Rosignoli, M.T., *et al.* Benzydamine protection in a mouse model of endotoxemia. *Inflamm. Res.* **46**(9), 332-335 (1997).
3. Alaiye, A., Kaya, E., Pınarbaşı, M.Ö., *et al.* An experimental comparison of the analgesic and anti-inflammatory effects of safflower oil, benzydamine HCl, and naproxen sodium. *J. Med. Food* **23**(8), 862-869 (2020).

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, 12/11/2023

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