

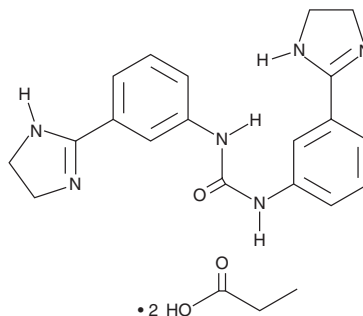
PRODUCT INFORMATION



Imidocarb (propionate)

Item No. 27053

CAS Registry No.: 55750-06-6
Formal Name: N,N'-bis[3-(4,5-dihydro-1H-imidazol-2-yl)phenyl]urea, dipropionate
MF: $C_{19}H_{20}N_6O \cdot 2C_3H_6O_2$
FW: 496.6
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 220, 260 nm
Supplied as: A crystalline solid
Storage: 4°C
Stability: ≥ 4 years



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

Laboratory Procedures

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of imidocarb (propionate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of imidocarb (propionate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Imidocarb is an antiprotozoal agent.¹ It is active against the protozoan parasite *B. bovis* ($IC_{50} = 0.87 \mu\text{g/ml}$).² Imidocarb (1.2 mg/kg) eliminates parasites in blood in a splenectomized lamb model of *B. ovis* infection and increases survival when administered at a dose of 2.4 mg/kg prior to infection.³ Formulations containing imidocarb have been used in the veterinary treatment of babesiosis and other protozoal infections.

References

1. Mosqueda, J., Olvera-Ramírez, A., Aguilar-Tipacamú, G., *et al.* Current advances in detection and treatment of babesiosis. *Curr. Med. Chem.* **19**(10), 1504-1518 (2012).
2. Rodriguez, R.I. and Trees, A.J. In vitro responsiveness of *Babesia bovis* to imidocarb dipropionate and the selection of a drug-adapted line. *Vet. Parasitol.* **62**(1-2), 35-41 (1996).
3. Sevinc, F., Turgut, K., Sevinc, M., *et al.* Therapeutic and prophylactic efficacy of imidocarb dipropionate on experimental *Babesia ovis* infection of lambs. *Vet. Parasitol.* **149**(1-2), 65-71 (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, 11/07/2022

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