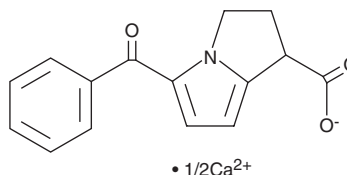


PRODUCT INFORMATION

Ketorolac (calcium salt)

Item No. 9003475

CAS Registry No.: 167105-81-9
Formal Name: 5-benzoyl-2,3-dihydro-1H-pyrrolizine-1-carboxylic acid, hemicalcium salt
MF: $C_{15}H_{12}NO_3 \bullet 1/2Ca$
FW: 274.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 246, 315 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Laboratory Procedures

Ketorolac (calcium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the ketorolac (calcium salt) in the solvent of choice, which should be purged with an inert gas. Ketorolac (calcium salt) is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Ketorolac (calcium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ketorolac (calcium salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ketorolac (calcium salt) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ketorolac is a non-steroidal anti-inflammatory drug (NSAID) and a non-selective COX inhibitor ($IC_{50} = 20$ nM for both COX-1 and COX-2).¹ It prevents increases in paw swelling, increases paw withdrawal latency in a hot-plate test, and decreases prostaglandin E_2 (PGE_2) levels in paw tissue in a mouse model of carrageenan-induced inflammation when administered at a dose of 30 mg/kg. Ketorolac is a racemic mixture containing the active (S)-ketorolac (Item No. 11348) and inactive (R)-ketorolac enantiomers. Formulations containing ketorolac have been used to manage postoperative pain and as ophthalmic solutions to treat ocular pain and inflammation.

Reference

1. Zhang, Y., Shaffer, A., Portanova, J., *et al.* Inhibition of cyclooxygenase-2 rapidly reverses inflammatory hyperalgesia and prostaglandin E_2 production. *J. Pharmacol. Exp. Ther.* **283**(3), 1069-1075 (1997).

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, 06/30/2020

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