

PRODUCT DATA SHEET



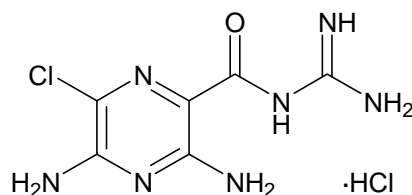
PRODUCT: Amiloride-HCl
3,5-Diamino-N-(aminoiminomethyl)-6-chloropyrimidinamide

CAS NO: 2016-88-8

CATALOG NO.: CA-200

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: Off-white solid (m.p. 288-293°C)

MOLECULAR FORMULA: C₆H₈ClN₇O·HCl MOLECULAR WEIGHT: 266.1

PURITY: 99% (TLC)

SOLUBILITY: Soluble in DMSO (25 mg/mL) and water (10 mM, warm)

STORAGE: Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.

APPLICATION NOTES: Selectively blocks low threshold (T-type) calcium channel ($K_D = 30 \mu\text{M}$) with no effect on L-type channels or tetrodotoxin Na⁺ channels.¹ Inhibits the epithelial tetrodotoxin insensitive Na⁺ channel ($IC_{50} = 1 \mu\text{M}$).²⁻⁴ Clinically useful K⁺ sparing diuretic.⁵

- REFERENCES:
1. C.-M.Tang *et al. Science* 1988 **240** 213
 2. S.Saribon-Sohraby and D.Benos *Am.J.Physiol.* 1986 **250** C175
 3. T.R.Kleyman and E.J.Cragoe, Jr. *J.Membrane Biol.* 1988 **105** 1
 4. T.Matainaho *et al. J.Endocrinol.* 1989 **123** 93
 5. H.L.Macfie *et al. Drug Intell.Clin.Pharmacol.* 1981 **15** 94

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.