

PRODUCT DATA SHEET



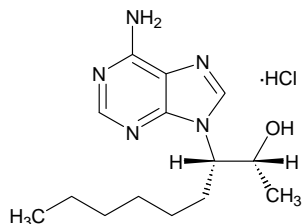
PRODUCT: EHNA
erythro-9-(2-Hydroxy-3-nonyl)adenine·HCl

CAS NO: 51350-19-7

CATALOG NO.: PD-129

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 183-184°C dec.)

MOLECULAR FORMULA: C₁₄H₂₃N₅O·HCl

MOLECULAR WEIGHT: 313.8

PURITY: 98%

SOLUBILITY: Soluble in water (25 mg/ml)

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

APPLICATION NOTES: Phosphodiesterase PDE II inhibitor (IC₅₀ = 0.8 μM).^{1,2} Also inhibits adenosine deaminase (K_i = 4 nM).³

REFERENCES: 1. T.Podzuweit *et al. Cell.Signal.* 1995 **7** 733
2. P.F.Mery *et al. Mol.Pharmacol.* 1995 **48** 121
3. M.Bessodes *et al. Biochem.Pharmacol.* 1982 **31** 879

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.