

# PRODUCT DATA SHEET



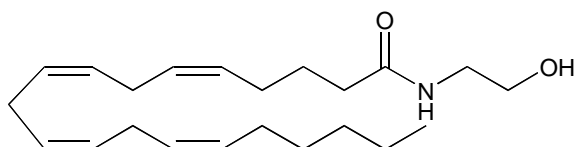
PRODUCT: Anandamide

CAS NO: 94421-68-8

CATALOG NO.: FA-017

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: Colorless oil

MOLECULAR FORMULA:  $C_{22}H_{37}NO_2$

MOLECULAR WEIGHT: 347.6

PURITY: 98% (TLC)

STORAGE: Store, as supplied, under an inert atmosphere at  $-80^{\circ}C$  for up to 1 year.

APPLICATION NOTES: After opening, protect from oxidation. Endogenous cannabinoid ligand.<sup>1</sup> Inhibits the binding of [ $^3H$ ]HU-243 to synaptosomal membranes ( $K_i = 52$  nM).<sup>2</sup> Binds to brain<sup>2</sup> and peripheral<sup>3</sup> cannabinoid receptors. Cannabimimetic activity.<sup>4,5</sup>

- REFERENCES:
1. W.A.Devane *Trends Pharmacol.Sci.* 1994 **15** 40
  2. W.A.Devane *et al. Science* 1992 **258** 1946
  3. S.Munro *et al. Nature* 1993 **365** 61
  4. E.Fride *et al. Eur.J.Pharmacol.* 1993 **231** 313
  5. J.N.Crawley *et al. Pharmacol.Biochem.Behav.* 1993 **46** 967

**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.**