

PRODUCT DATA SHEET



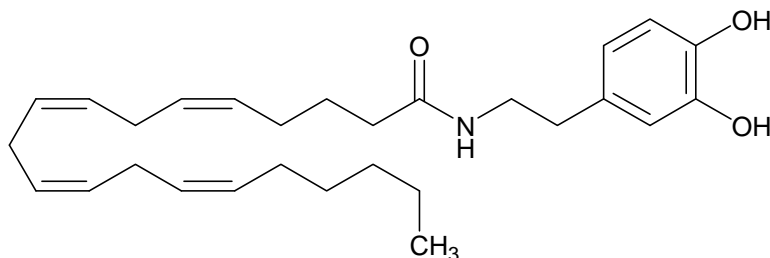
PRODUCT: N-Arachidonoyldopamine
NADA

CAS NO: 199875-69-9

CATALOG NO.: VR-104

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: Viscous oil

MOLECULAR FORMULA: $C_{28}H_{41}NO_3$

MOLECULAR WEIGHT: 439.6

PURITY: 98% (TLC: 5% Methanol/methylene chloride; $R_f = 0.20$)

SOLUBILITY: Soluble in DMSO

STORAGE: Store, as supplied, at $-20^{\circ}C$ for up to 1 year. Store solutions at $-20^{\circ}C$ for up to 3 months.

APPLICATION NOTES: Endogenous ligand for the VR1 receptor ($EC_{50} = 63$ nM).^{1,2} Selective agonist at cannabinoid CB_1 versus CB_2 receptors ($K_i = 0.25$ and 12 μM respectively).³ Product is subject to oxidation. Protect from air. Store under an inert atmosphere.

- REFERENCES:
1. C.J.Chu *et al. J.Biol.Chem.* 2003 **278** 13633
 2. S.M.Huang *et al. Proc.Natl.Acad.Sci.USA* 2002 **99** 8400
 3. T.Bisogno *et al. Biochem.* 2000 **351** 817

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.