

# PRODUCT DATA SHEET



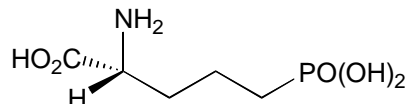
PRODUCT: D(-)-2-Amino-5-phosphonovaleric acid  
D-AP5

CAS NO: 79055-68-8

CATALOG NO.: EA-104

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 233°C)

MOLECULAR FORMULA: C<sub>5</sub>H<sub>12</sub>NO<sub>5</sub>P

MOLECULAR WEIGHT: 197.1

PURITY: 99%

SOLUBILITY: Soluble in water

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

APPLICATION NOTES: Potent NMDA antagonist. Active form of AP5.<sup>1</sup>

REFERENCES: 1. J.C.Watkins *et al. Ann.Rev.Pharmacol.Toxicol.* 1981 **21** 165  
2. J.J.Hansen and P.Krogsgaard-Larson *Med.Res.Rev.* 1990 **10** 55

**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.**