

PRODUCT DATA SHEET



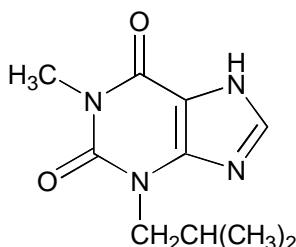
PRODUCT: IBMX
3-Isobutyl-1-methylxanthine

CAS NO: 28822-58-4

CATALOG NO.: PD-140

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 199-201°C)

MOLECULAR FORMULA: C₁₀H₁₄N₄O₂

MOLECULAR WEIGHT: 222.2

PURITY: 99% (TLC)

SOLUBILITY: Soluble in DMSO and ethanol

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

APPLICATION NOTES: A non-specific inhibitor of phosphodiesterases (IC₅₀ = 2-50 μM).¹ Also possesses adenosine agonist activity.^{2,3}

- REFERENCES:
1. J.A. Beavo *et al. Mol. Pharmacol.* 1970 **6** 597
 2. Bruns *et al. Mol. Pharmacol.* 1986 **29** 331
 3. Coffin *et al. Eur. J. Pharmacol.* 1989 **170** 35

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.