

Milrinone

Catalog #: 27216

Lot #: 120608

Size: 10 mg

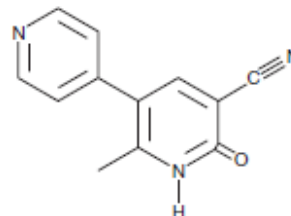
Structure:

CAS Registry #: 78415-72-2

Purity: ≥98%

Chemical Formula: C₁₂H₉N₃O

Molecular Weight: 211.2



Description: Milrinone, also known as WIN 47203 and Primacor[®], is an inhibitor of PDE3. It is less effective against other PDE isoforms. Milrinone has positive inotropic and vasodilatory effects *in vivo*.

Appearance: A crystalline solid

Solubility: Soluble in DMSO and DMF, purged with an inert gas. Solubility in DMSO and DMF is ~0.3 mg/ml. Do not store aqueous solutions for more than one day.

Biological Activity: Milrinone inhibits PDE3A and PDE3B with IC₅₀ values of 0.45 μM and 1.0 μM, respectively.

Storage/Stability: Store at or below -20°C for up to two years.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Sudo, T., *et al.*, *Biochem Pharmacol.* 2000;**59**:347-56.
2. Lakshminrusimha, S., *et al.*, *Pediatr. Crit. Care Med.* 2009;**10**(1):106-12.
3. Simkova, V., *et al.*, *Crit. Care.* 2007;**11**(3):139.
4. Maskin, C.S., *et al.*, *Circ. J.* 1983;**67**:1065-70.