

Instruction Manual

| | |
|---|---|
| Catalog Number | PK-CA577-1041-1G |
| Description | A very active antibiotic against many yeasts and fungi. Inhibits protein synthesis in cytoplasmic eukaryotes but not in prokaryotes. Completely inhibits many molds but is noninhibitory at 100 µg/ml to many species of pathogenic bacteria. It inhibits chain initiation as well as chain elongation by acting on the 60S subunit of the eukaryote ribosome, interacting directly with enzyme translocase. Cycloheximide is used as an inhibitor to study cell-free protein biosynthesis in eukaryotes and also used to block ribosome-dependent in vivo polypeptide synthesis. It is an inducer of apoptosis in many cell lines (1). |
| Quantity | 1 g |
| Sequence / Molecular Weight / Molecular Formula | 281.3 Da; C ₁₅ H ₂₃ NO ₄ |
| Purity | >95% |
| Appearance / Formulation / Solubility | Solid |
| Storage & Stability | Store powder at desiccated at -20°C for up to 1 year. Store solution at -20°C for up to 3 months. Rapidly decomposed in alkaline conditions. |
| Applications | see Description |
| References | 1. Grand, R.J., et al. (1995) Exp. Cell Res. 218:439. 2. Lotem, J. and L. Sachs. (1995) Cell Growth Diff. 6:647. 3. Chow, S.C., et al. (1995) Exp. Cell Res. 216:149. 4. Alten, J., et al. (1995) Cell Immunol. 161:98. |
| Caution | NA |

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.
