

Product Information

| | |
|------------------------------|--|
| Description: | The NP-40 Cell Lysis Buffer is intended for the lysis of mammalian adherent and suspension-cultured cells. This buffer is suitable for extraction of cytoplasmic proteins in mild, non-denaturing conditions. |
| Formulation: | 50 mM Tris HCl, pH 7.4, 150 mM NaCl, 1 mM EDTA, 1 mM EGTA, and 1% IGEPAL CA-630 |
| Quality: | Sterile filtered (0.2 μ m) |
| Storage/Stability: | Upon receipt, store at 4°C. Stable for 6 months from date of receipt, when stored as directed. |
| Instructions for Use: | This is a ready-to-use buffer. Mix well before use. This buffer does not contain protease or phosphatase inhibitors. If desired, add appropriate inhibitors (such as protease inhibitors, phosphatase inhibitors, ADP-ribosylation inhibitors etc.) immediately before use to prevent protein proteolysis or to maintain phosphorylation and/or ribosylation status of proteins. |
| Shipping Temperature: | 4°C (gel packs) |