

Jurkat cytoplasmic extract

Catalog No: 36402

Contents: 200 µg of Jurkat cytoplasmic extract at 5.0 µg/µl

Composition:

Cytoplasmic extracts are supplied in dilution buffer (10mM Hepes (pH 7.9), 10 mM KCl, 0.1mM EDTA, 0.1mM EGTA, 0.5 mM PMSF, 1 mM DTT, 1ug/ml Aprotinin and 1ug/ml Leupeptin).

Quality Control:

Extracts have been quality control tested by Western blot .

Storage:

To ensure stability, extracts should be stored at -80°C.

We recommend aliquoting the extracts into single-use fractions and then storing them at -80°C. This eliminates repeated freeze/thaw cycles.

Guarantee:

This product is guaranteed for 6 months from date of arrival.