

Nb2 nuclear extract

Catalog No: 36105

Contents: 2 x 100 µg of Nb2 nuclear extract at 2.5 µg/µl

Composition:

The Nb2 nuclear extract derived from rat lymphocytes have been collected in Lysis Buffer. Lysis Buffer consists of 20 mM Hepes pH 7.9, 100 mM KCl, 20% glycerol, 1 mM MgCl₂, 0.5 mM DTT and 0.5 mM PMSF. The protein content has been determined by a Bradford-based assay.

Quality Control:

Each lot has been tested for STAT activation by using TransAM™ STAT Family Kits.

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.