

F9 nuclear extract (Retinoic acid treated)

Catalog No: 36113

Contents: 2 x 100 µg of F9 nuclear extract (Retinoic acid treated) at 2.5 µg/µl

Composition:

Nuclear extracts are supplied in dilution buffer (20mM Hepes (pH 7.9), 100 mM KCl, 1mM MgCl₂, 20% glycerol, 0.5 mM PMSF and 0.5 mM DTT). Cells were treated with 1 µM Retinoic acid for 48 hours. Retinoic acid is refreshed after 24 hours. Cells must be grown on plates coated with 0.1% gelatin.

Quality Control:

Each lot has been tested for GATA-4 activation by using TransAM™ 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.