
MCF-7 Nuclear Extract

(Catalog No. X12002)

Concentration

Supplied in four vials, each containing 250 µg nuclear extract in 50 µl buffer

Size

1000 µg

Storage Buffer

Provided in 20 mM HEPES (pH 7.9), 20% v/v glycerol, 0.1 M KCl, 0.2 mM EDTA, 0.5 mM PMSF and 0.5 mM DTT

Storage

Extracts should be stored at -80°C and repeated freezing and thawing should be avoided.

Extract Origin

Human nuclear extract; mammary adenocarcinoma cells

Applications

Suitable for use in Gel Shift and Western Blotting assays

Ordering Information:

Products	Size	Cat. No.
MCF-7 Nuclear Extract	1000 µg	X12002 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.