

---

## DU-145 Nuclear Extract

(Catalog No. X12000)

---

### 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; prostate carcinoma cells

### Applications

Suitable for use in Gel Shift and Western Blotting assays

### Ordering Information:

Products	Size	Cat. No.
DU-145 Nuclear Extract	1000 µg	X12000 -1

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