

PRODUCT NAME**Unblocked protein A beads**

Cat. No: kch-503-880	Format: 880 μ l
Cat. No: kch-503-028	Format: 2.8 ml
Cat. No: kch-503-008	Format: 8 ml

Description

The unblocked protein A beads are agarose beads that have been extensively validated in chromatin immunoprecipitation assays (ChIP). These beads are intended for isolation of immune complexes (chromatin and specific antibody) in ChIP experiments. The beads are suitable for immunoprecipitation of rabbit polyclonal Abs, mouse IgG2a, IgG2b and IgA, guinea pig IgG, dog IgG, pig IgG. The beads should be washed before use.

Format

Supplied as a suspension 1:4 in PBS with 20% of ethanol. Keep the beads in liquid suspension during storage as drying will result in reduced performance.

Storage

Store at 4°C. Do not freeze. Stable for 1 year.

Precautions

This product is for research use only. Not for use in diagnostic or therapeutic procedures.