

NB500-213 Protocol

Immunoprecipitation protocol specific for Smac Antibody (NB500-213)

SMAC/Diablo Antibody (78-1-118): https://www.novusbio.com/products/smac-diablo-antibody-78-1-118_nb500-213
Procedure Guide for NB 500-213 Monoclonal anti-Survivin (clone 58.22)

Immunoprecipitation Procedure

1) Lyse cells plated in a 60mm dish:

a) 300 ul CHAPS buffer [50mM Tris-HCl, pH 7.5/50mM NaCl/1mM EDTA/1% NP-40/0.1% CHAPS/1mM NaVO₄/1mM PMSF]

b) rock for 20 minutes at 4C

2) Harvest lysate and spin down the insoluble material at 14K rpm.

3) Collect soluble fraction.

4) Pre-clear lysate with 40 ul of 50:50 slurry of Protein A beads, rocking for 1 hour at 4C.

5) Spin down beads at 2K rpm, at 4C.

6) Collect pre-cleared lysate.

7) Incubate lysate with 5-7ug of anti-Survivin (NB 500-213) overnight, rocking at 4C.

8) Add 50 ul of Protein A 50:50 slurry for 2 hours, rocking at 4C.

9) Wash beads with 200 ul of CHAPS buffer, three times.

10) Denature immune complex by adding 2x Sample Buffer, containing 2-ME.

11) Boil for 10 minutes and load onto an SDS-gel.