

## DIMENSIONS OF TOPSEAL-A PLUS (PN 6050185)

### Material

Clear polyester film coated with a clear pressure sensitive acrylic adhesive specially designed for the rapid and easy sealing of polystyrene microplates. The perforated end tabs (see picture) allow easy handling, accurate positioning and quick removal of the tabs after sealing. Each box contains 100 seals and an instruction card.

### Applications

TopSeal-A Plus can be used with microplates for radioisotopic and luminescence analysis on the TopCount or MicroBeta counters from PerkinElmer.

### Dimensions

Length of seal with tabs	152	mm
Length of seal without tabs	120	mm
Width of seal	78	mm
Seal thickness without backing	0.07	mm

