

Research use only. Not for use in diagnostic procedures.

## Dimensions of Yellow DELFIA Plate

<b>Part numbers:</b>	PN AAAND-0001	Transparent yellow, high protein binding affinity (HB)
	PN AAAND-0004	Transparent yellow, anti-rabbit IgG coated
	PN AAAND-0005	Transparent yellow, streptavidin-coated

**Material and treatment or coating:**

Transparent yellow polystyrene microplate with 96 flat-bottom wells. Plates have high protein binding affinity and are available uncoated, coated with streptavidin or coated with anti-rabbit IgG's. Plates have a transparent yellow color by added UV absorber for reduced background in DELFIA assays.

**Packaging:**

AAAND-0001 is a case with 60 microplates packed as 12 stacks of 5 plates each. AAAND-0004 and AAAND-0005 is a box with 10 individual packaged coated plates.

**Dimensions --- Customer Drawing:**

