

## Product Information

### 6X DNA Loading Buffer

Catalog Number	Packaging Size
D016	5 X 1 mL

#### Storage upon receipt:

- Room temperature

### Product Description

**DNA Loading Buffer** is a 6X solution of 30mM EDTA, 36% (v/v) Glycerol, 0.05% (w/v) Xylene Cyanol FF, 0.05% (w/v) Bromophenol Blue for use in agarose gel electrophoresis of nucleic acids.

### Application

1. Mix 5 volumes of DNA sample with 1 volume of **DNA Loading Buffer (6X)**. Vortex briefly.
2. Load on the gel, and run electrophoresis.

### Related Products

Catalog No.	Product
D012	<b>SafeGreen™ Loading Dye with DNA Loading Buffer (6X)</b>
D013	<b>SafeRed™ Loading Dye with DNA Loading Buffer (6X)</b>