



## Blue Gel Loading Buffer

Loading buffer for agarose or polyacrylamide gels

Cat. No.	Amount
PCR-254-bl	5 x 1,8 ml

**For in vitro use only!**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at 4 °C

**Shelf Life:** 12 months

**Form:** Liquid

**Concentration:** 6x conc.

### Applications:

Add one part of Gel Loading Buffer to 5 parts of the DNA sample as shown in the table below. Spin down the tubes and vortex gently to achieve homogeneity.

DNA sample	5 µl	10 µl	20 µl	50 µl
6x Gel Loading Buffer	1 µl	2 µl	4 µl	10 µl

### Description:

Jena Bioscience Gel Loading Buffers are formulated to facilitate loading of DNA containing samples into the wells of agarose and polyacrylamide gels. The buffers contain tracking dyes as indicator for DNA fragment migration. In addition, they contain glycerol to add density and EDTA to inhibit nuclease activities.

The buffers are optimized for loading of DNA fragments in a size range of:

Cat. No.	Gel Loading Buffer	fragment size
PCR-254-bl	Blue Gel Loading Buffer	larger than 500 bp
PCR-254-gr	Green Gel Loading Buffer	from 100 to 2000 bp
PCR-254-or	Orange Gel Loading Buffer	smaller than 500 bp

### 6x Gel Loading Buffer:

60 mM Tris-HCl (pH 7.5), 60 mM EDTA, 50 % (w/v) Glycerol, bromophenol blue (blue buffer), xylene cyanol FF (blue and green buffer), orange G (orange buffer) and tartrazine (green buffer)

DNA sample	5 µl	10 µl	20 µl	50 µl
6x Gel Loading Buffer	1 µl	2 µl	4 µl	10 µl



## Blue Gel Loading Buffer

Loading buffer for agarose or polyacrylamide gels

### Fragment Separation on Agarose Gels:

DNA fragment size	agarose gel conc.	xylene cyanol FF running at approx.	bromo-phenol blue running at approx.	orange G running at approx.
20-400 bp	3.6 %	280 bp	40 bp	2 bp
50-1000 bp	3.0 %	500 bp	60 bp	2 bp
100-2000 bp	2.4 %	900 bp	100 bp	3 bp
200-4000 bp	1.8 %	1800 bp	40 bp	5 bp
0.5-10 kb	1.2 %	4.5 kb	0.5 kb	10 bp
1-30 kb	0.6 %	12 kb	1.2 kb	100 bp