



175PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# Gel Loading Dye

## For Loading DNA & RNA Samples

(Cat. # 786-100 to 786-107)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

Gel Loading Dyes are ready to use dyes for running agarose gel electrophoresis of DNA and RNA. The dyes for DNA are Ficoll based and are available as Glow™ Loading Dye and Universal Loading Dye. There is no need to add ethidium bromide to the running buffer as Glow™ Loading Dyes contain ethidium bromide, which reduces exposure and many of the problems associated with it. Glow™ Loading Dyes provide intense DNA bands with little background or band distortion and can be used with any type of agarose or acrylamide gels. Also available are Universal Loading Dyes with an option of added SDS.

RNA Loading Dyes are formamide based and are supplied as 2X concentration. The Glow™ RNA Dye is ideal for quick screenings of RNA preps. There is no need to use toxic formaldehyde agarose gels for checking RNA integrity with Glow™ RNA Dye.

## ITEM(S) SUPPLIED

The following tables show the general composition of the various loading dyes.

### Universal Dyes [6X]

Cat. #	Description	Bromophenol Blue	Xylene Cynol	SDS	Size
786-100	BromoBlue Universal Dye	Yes	-	-	1.5ml
786-101	CyanoBlue Universal Dye	Yes	Yes	-	1.5ml
786-102	CleanAway™ Universal Dye	Yes	Yes	Yes	1.5ml

### Glow™ Dyes [6X]

Cat. #	Description	Bromophenol Blue	Xylene Cynol	SDS	Ethidium Bromide	Size
786-103	Glow™ BromoBlue Dye	Yes	No	No	Yes	1.5ml
786-104	Glow™ CyanoBlue Dye	Yes	Yes	No	Yes	1.5ml
786-105	Glow™ CleanAway™ Dye	Yes	Yes	Yes	Yes	1.5ml

### RNA Loading Dyes [2X]

Cat. #	Description	Bromophenol Blue	Xylene Cynol	Ethidium Bromide	Size
786-106	Universal RNA Dye	Yes	Yes	No	1.5ml
786-107	Glow™ RNA Dye	Yes	Yes	Yes	1.5ml

## STORAGE

It is shipped at ambient temperature. Store Gel Loading Dye at 4°C, upon arrival.

## PROTOCOL

### **6X Universal & Glow Dyes**

Dilute one part 6X Dye solution into five parts of sample solution to give a final concentration of 1X Dye solution. The sample is then ready to load to a gel.

For Example: 10µl sample and 2µl 6X Dye Solution.

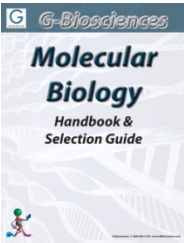
### **2X RNA Dyes**

Mix equal volumes of 2X Dye Solution and RNA sample solution to give a final concentration of 1X Dye solution. The sample is then ready to load on a gel.

For example: 10µl sample and 10µl 2X RNA Dye Solution.

## RELATED PRODUCTS

Download our Molecular Biology Handbook.



<http://info.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

Last saved: 8/1/2012 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)