

Gel Loading Dye

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**.

KIT COMPONENTS

1 Tube containing 1.5ml dye for each catalog number as follows:

Catalog #	Description	Bromophenol Blue	Xylene Cynol	SDS	Ethidium Bromide	Size
<u>6X Universal Dyes</u>						
786-100	BromoBlue Universal Dye	Yes	No	No	No	1.5ml x 1
786-101	CyanoBlue Universal Dye	Yes	Yes	No	No	1.5ml x 1
786-102	CleanAway™ Universal Dye	Yes	Yes	Yes	No	1.5ml x 1
<u>6X Glow™ Dyes</u>						
786-103	Glow™ BromoBlue Dye	Yes	No	No	Yes	1.5ml x 1
786-104	Glow™ CyanoBlue Dye	Yes	Yes	No	Yes	1.5ml x 1
786-105	GlowCleanAway™ Dye	Yes	Yes	Yes	Yes	1.5ml x 1
<u>2X RNA Dyes</u>						
786-106	Universal RNA Dye	Yes	Yes	No	No	1.5ml x 1
786-107	Glow™ RNA Dye	Yes	Yes	No	Yes	1.5ml x 1

STORAGE

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 + 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 + 10µl 2X RNA Dye Solution.

RELATED PRODUCTS

1. **femtoLUCENT™ -A Chemiluminescence Detection System (Cat # 786-10)** - Genotech has developed the femtoLUCENT kit, which is based on isoluminol-HRP-chemiluminescence reaction. The chemiluminescence light emission can be recorded by a short exposure to blue-light sensitive autoradiography film. femtoLUCENT™ allows detection of sub-pico gram quantity of antigens with lower background.
2. **Ready-to-Screen Tissue Blots for Western Analysis:** G-Biosciences has put together pre-made Ready-to-Screen Tissue Blots of total proteins extracted from hard-to-obtain tissues from human (normal, tumor and region specific), mouse, rats and cell lines etc.
3. **Western Re-Probe™ (Cat # 786-119)** - Western Re-Probe (5X) kit provides a buffer for stripping and re-probing western blot membranes.
4. **NAP-Blocker™ (Cat # 786-190)** – G-Biosciences's NAP-Blocker™ (supplied as 2X) is a blocking agent containing non-animal proteins for improved assay sensitivity, clean background and a high signal to background ratio. Simple to use and better results than milk powder preparation as it ensures uniform blocking without non-specific binding.

For more information on these and other products, please visit our website at www.GBiosciences.com