

# TREVIGEN® Product Data

*For Research Use Only. Not For Use In Diagnostic Procedures*

---

## 5X Orange G Loading Buffer

Cat# 9850-250-6

| Contents                   | Catalog#   | Size       | Storage    |
|----------------------------|------------|------------|------------|
| 5X Orange G Loading Buffer | 9850-250-6 | 250 µl x 6 | Room Temp. |

**Description:** Trevigen's **5X Orange G Loading Buffer** is an ideal sample loading buffer for PCR assays and DNA electrophoresis on agarose gels. The high density of the loading buffer ensures that the DNA samples load easily and remain at the bottom of the preformed wells of agarose gels. The orange colored dye front migrates at the 10 - 20 base pair range and does not interfere with the migration of DNA fragments. Thus, any obscuring of smaller fragments is prevented. Additionally, 5X Orange G Loading Buffer does not interfere with viewing DNA fragments under ultraviolet light in ethidium bromide stained gels.

Trevigen's Orange G Loading Buffer is provided at a 5X concentration and does not require refrigeration. It is also recommended for use with TreviGel™ 500 and TreviGel™ 5000, and for analysis of small DNA products of PCR assays.

©2007 Trevigen, Inc. All rights reserved. Trevigen is a registered trademark of Trevigen, Inc.

---

# TREVIGEN®

8405 Helgerman Court, Gaithersburg, MD 20877 USA  
 Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800  
 Fax: 301-560-4973 • e-mail: info@trevigen.com • www.trevigen.com

**5X Orange G  
Loading Buffer**

Catalog #:9850-250-6

Storage: Room Temperature

**TREVIGEN®**

1-800-873-8443