## Detection of the beta-galactosidase Reporter Gene in Transfected Eukaryotic Cells

This protocol is for the Detection of the beta-galactosidease Reporter Gene in Transfected Eukaryotic Cells

**10X PBS buffer (pH 7.4):** 1.37 M NaCl, 0.27 M KCl, 1 M Na, HPO<sub>4</sub>, 0.02 M K, HPO<sub>4</sub>.

Fixation buffer (pH 7.4): 1X PBS buffer and 0.25% glutardialdehyde.

**Staining buffer,** prepare immediately before use as follows:

Stock solutions	Volume per 10 ml staining buffer	Final concentration
1 M MgCl2	20 μΙ	2 mM
0.5 M K4Fe(CN)63H2O	100 μΙ	5 mM
0.5 M K3Fe(CN)6	100 μΙ	5 mM
X-Gal Solution, ready-to-use	500 µІ	1 mg/ml
10X PBS buffer (pH 7.4)	9.28 ml	diluted 10-fold

## Staining procedure:

- 1. Wash the cells twice with cold 1X PBS buffer. Adhered cells can be washed in the transfection plates, suspension cells should be pelleted before washing.
- Fix the cells with Fixation buffer for 10 minutes at room temperature while gently rocking the plate.
   Use 150 μl of the Fixation buffer for each well of a 24-well plate.
- 3. Wash the cells twice with cold 1X PBS buffer.
- 4. Stain the cells with freshly prepared Staining buffer for 2-20 hours at 37°C. Use 200 μl of Staining buffer for each well of a 24-well plate.
- 5. Count dark blue cells.

## thermoscientific.com

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

North America
Technical Services:
techservice.genomics@thermofisher.com
Customer Services:
customerservice.genomics@
thermofisher.com
Tel 800 235 9880
Fax 800 292 6088

Technical Services: techservice.emea.genomics@ thermofisher.com Customer Services: customerservice.emea.genomics@ thermofisher.com

