

CyQUANT™ Direct Cell Proliferation Assay

Catalog No. C35011, C35012

Pub. No. MAN0018318

Rev. A.0

Upon receipt, kit can be stored at room temperature for 6 months.

1. For each microplate, prepare 12 mL of 2X Detection Reagent as follows:

PBS, HBSS, or cell culture media (\pm serum)	11.7 mL
CyQUANT™ Direct Nucleic Acid Stain	48 μ L
CyQUANT™ Direct Background Suppressor	240 μ L
Total Volume	12 mL

2. Add an equal volume of 2X Detection Reagent to cells in cell culture media (\pm serum).
3. Incubate the cells with the Detection Reagent for 60 minutes at 37°C.
4. Read the fluorescence from bottom of the well with standard green filter sets or wavelengths (ex/em maxima 508/527 nm).

For Research Use Only. Not for use in diagnostic procedures.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product you accept the terms and conditions of all applicable Limited Use Label Licenses.

Manufacturer: Life Technologies Corporation | 29851 Willow Creek Road | Eugene, OR 97402

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit thermofisher.com/support
thermofisher.com

ThermoFisher
 S C I E N T I F I C