

Product Number
829-07921

Quantity: 25 µL

Storage: -20°C

Revised: July 2011

Updates available at:
<http://biosupport.licor.com>

**Limitation of Liability and
Limited Use Label License**

LI-COR IRDye® Infrared Dyes and reagent products are offered for research purposes only and are not intended for human therapeutic or diagnostic use. The purchase of this product conveys to the buyer the non-transferable right to use the amount of product purchased and the components of the product in research conducted by the buyer (whether the buyer is a not-for-profit, academic or for-profit entity). The buyer shall not sell or otherwise transfer this product, its components, or materials made there from to any third party. Buyer shall not use this product or its components for commercial purposes. The term "commercial purposes" shall mean any activity by a party for consideration and may include, but is not limited to, use of the product or its components (i) in manufacturing, (ii) to provide a service, information or data, (iii) for therapeutic, diagnostic or prophylactic purposes, or (iv) for resale, whether or not such product or its components are resold for use in research. The use of this product by the buyer constitutes agreement with the terms of this limited use label license for LI-COR IRDye infrared dyes and reagent products. Inquiries regarding the licensing of one or more IRDye infrared dyes should be submitted by e-mail to busdev@licor.com.

LI-COR BIOSCIENCES DOES NOT PROVIDE RESEARCH ADVICE OR DETERMINE OR RECOMMEND ANY POTENTIAL USES FOR IRDYE INFRARED DYES AND REAGENT PRODUCTS. LI-COR BIOSCIENCES MAKES NO WARRANTIES OF ANY KIND, EITHER EXPRESS OR IMPLIED, AS TO ANY MATTER INCLUDING, BUT NOT LIMITED TO, WARRANTY OF FITNESS FOR PURPOSE, OR MERCHANTABILITY OR RESULTS OBTAINED FROM USE OF IRDYE INFRARED DYES. IN NO EVENT SHALL LI-COR BE LIABLE FOR LOST PROFITS, CONSEQUENTIAL, EXEMPLARY, SPECIAL, DIRECT, INCIDENTAL, OR PUNITIVE DAMAGES, OR ATTORNEY FEES, EVEN IF LI-COR HAD BEEN ADVISED OF, KNEW, OR SHOULD HAVE KNOWN, OF THE POSSIBILITIES THEREOF. NO EMPLOYEE, AGENT OR REPRESENTATIVE OF LI-COR HAS THE AUTHORITY TO BIND LI-COR TO ANY ORAL REPRESENTATION OR WARRANTY EXCEPT AS SPECIFICALLY SET FORTH HEREIN.

© 2011 LI-COR, Inc. LI-COR is an ISO 9001 registered company. The Odyssey Imager, Aeriis Imager, and the IRDye reagents are covered by U.S. and foreign patents and patents pending. LI-COR, IRDye, Aeriis, and Odyssey are registered trademarks of LI-COR, Inc. All other trademarks belong to their respective owners.

Doc #988-12279

LI-COR

Biosciences

4647 Superior Street • P.O. Box 4000
Lincoln, Nebraska 68504 USA
North America: 800-645-4267
International: 402-467-0700
FAX: 402-467-0819

LI-COR GmbH Germany, Serving Europe,
Middle East and Africa: +49 (0) 6172 17 17 771
LI-COR UK Ltd. UK, Serving UK, Ireland, and
Scandinavia: +44 (0) 1223 422104
All other countries, contact LI-COR Biosciences
or a local LI-COR distributor:
<http://www.licor.com/distributors>

www.licor.com

IRDye® 700 p53 Consensus Oligonucleotide

p53 Consensus Oligonucleotide¹

5' -- TAC AGA ACA TGT CTA AGC ATG CTG GGG ACT --3'
3' -- ATG TCT TGT ACA GAT TCG TAC GAC CCC TGA --5'

* Underlined nucleotides are the binding site

Introduction

Gel shift assays or electrophoretic mobility shift assays (EMSA) provide a simple method to study DNA-protein interactions. This assay is based on the principle that a DNA-protein complex will have a different mobility during electrophoresis than unbound DNA. These shifts can be visualized on a native acrylamide gel using labeled DNA to form the DNA-protein binding complex. The Aeriis® and Odyssey® family of imaging systems offer a quick and easily adapted alternative method to radioisotopic and chemiluminescent detection methods for EMSA analysis and visualization.^{2,3}

A DNA oligonucleotide end-labeled with IRDye 700 is a good substrate for protein binding. IRDye DNA detection is linear within a 50-fold dilution range from 9.1 fmol to 0.18 fmol. Additional benefits include no hazardous radioisotopes, no gel transfer to membrane or gel drying, no chemiluminescent substrate reagents, and no film exposure. Following electrophoresis, the gel can be imaged while in the glass plates. If necessary, the gel can be placed back in the electrophoresis unit and run longer.

Existing mobility shift assay protocols can be easily transformed into infrared assays by replacing the existing DNA oligonucleotides with IRDye oligonucleotides. The binding conditions and electrophoresis conditions will remain the same as with any other EMSA detection method.

Electrophoretic Mobility Shift Assay

A universal binding condition that applies to every protein-DNA interaction is not recommended, since binding conditions will be specific for each protein-DNA interaction. Thus, the user should establish the conditions of the binding reaction for each protein-DNA pair. Binding buffer should be the same for a specific DNA-protein complex as with any other mobility shift detection method used.

For IRDye 700 p53, the following binding reaction is a good starting point:

Reaction	µL
10X Binding Buffer (100 mM Tris, 500 mM KCl, 10 mM DTT; pH 7.5)	2
Poly (dI•dC) 1 µg/µL in 10 mM Tris, 1 mM EDTA; pH 7.5	1
25 mM DTT/2.5% Tween® 20	2
100 mM MgCl ₂	1
Water	12
IRDye 700 p53	1
HeLa 4 hour Serum Response nuclear extract (Positive control) (5 µg/µL)	1
Total	20

After the addition of the DNA to the protein-buffer mix, reactions are incubated to allow protein binding to DNA. A typical incubation condition is 20-30 minutes at room temperature. Since IRDye infrared dyes are somewhat sensitive to light, it is best to keep binding reactions in the dark during incubation periods (e.g., put tubes into a drawer or simply cover the rack containing tubes with aluminum foil). After the incubation period, Orange Loading Dye 10X (LI-COR, P/N 927-10100) is added to the binding reaction for electrophoresis.

IMPORTANT: It is critical **NOT** to use any blue loading dye (e.g., Bromophenol blue), as this will be visible on the Odyssey® image. It is highly recommended that Orange Loading Dye 10X (LI-COR, P/N 927-10100) be used instead.

NOTE: In some cases, we observed that DNA control reactions (no protein) have lower signal than the reactions containing the protein. This may be due to lower stability of the dye in certain buffer conditions. The addition of a final concentration of 2.5 mM DTT and 0.25% Tween® 20 to all reactions reduces this phenomenon.

Gel electrophoresis of the DNA-protein complex is done using a polyacrylamide gel composed of Tris-acetate, Tris-borate, or Tris-glycine-EDTA gel and buffer at 10 V/cm at room temperature or at 4°C in the dark (simply put a cardboard box over the electrophoresis apparatus).

Storage

Store at -20°C protected from light; stable for 6 months from date of shipment.

References

1. Kastan, M.B., Zhan, Q., El-Deiry, W.S., Carrier, F., Jacks, T., Walsh, W.V., Plunkett, B.S., Vogelstein, B., and Fornace, A.J.Jr.. 1992. *Cell* 71:587-597.
2. Li, Y., F. Ahmed, S. Ali, P. A. Philip, O. Kucuk, and F. H. Sarkar, 2005. *Cancer Res.* 65:6934-6942.
3. Geddie, M. L., T. L. O'Loughlin, K. K. Woods, and I. Matsumura, 2005. *J. Biol. Chem.* 280: 35641-35646.