



Product Announcement

The following reagent kits are for use with Agilent Gene Expression microarrays:

-
- | | |
|------------------|---|
| 5190-0442 | Quick Amp Labeling Kit, One-Color
Labeling reagents, including Cyanine 3-CTP; sufficient for 20 reactions, 20 arrays |
| 5190-0444 | Quick Amp Labeling Kit, Two-Color
Labeling reagents, including Cyanine 3- and Cyanine 5-CTP; sufficient for 20 reactions, 10 arrays |
| 5188-5282 | One color RNA Spike-in Kit
Reagents sufficient for up to 200 arrays |
| 5188-5279 | Two color RNA Spike-in Kit
Reagents sufficient for up to 200 arrays |
| 5188-5242 | Gene Expression Hybridization Kit with Hi-RPM Hyb Buffer
Reagents sufficient for 10 slides |
| 5190-0404 | Hi-RPM Gene Expression Hybridization Kit, Large Volume
Reagents sufficient for 100 slides |
| 5188-5327 | Gene Expression Wash Buffer Kit
Wash Buffers 1 & 2 sufficient for up to 80 slides |
| 5185-5979 | Stabilization & Drying Solution, 500 mL
Sufficient for up to 50 slides |
-

Products are for research use only. Not for use in diagnostic procedures.

Printed in USA

April 2008

Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610 USA



G4140-90080



Agilent Technologies