



STOP SOLUTION FOR GENOTYPING

Pack Size:	500 μ l	829-04459
	4 ml	830-05629
	200 ml	830-05630

Storage: -20 °C

RECOMMENDATIONS

- Aliquot the stop solution into several tubes of smaller volume.
- Store the tubes at -20 °C in the dark.
- Keep the stop solution on ice and return to the freezer after use.

DIRECTIONS FOR USE:

For standard 10 μ l reactions:

- After thermal cycling, add 5.0 μ l stop solution to each reaction.
- Denature the samples at 95 °C for 2-3 minutes and place on ice until ready to load the gel.

Note: The above volumes are guidelines. Please refer to LI-COR protocols in the genotyping manual for more information.

LI-COR[®]
Biotechnology Division

LI-COR, Inc. • 4308 Progressive Avenue • Lincoln, Nebraska 68504 USA
e-mail: biohelp@bio.licor.com • www.licor.com • Phone: 402-467-0700 • FAX: 402-467-0819
1-800-645-4267 (Toll free U.S. & Canada)

All brands/product names are trademarks of their respective owners.