

Additional cell vials

Terms and conditions

Version 24J02-NJ

DISCLAIMER

An “Additional vial” at the discounted price cannot be purchased alone and must be purchased simultaneously with a regular-priced vial of the same cell line.

CONTENTS

- **3-7 x 10⁶ cells** in a cryovial or shipping flask.

Note: If cells provided in a cryovial are not frozen upon arrival, contact InvivoGen immediately.

Important: Additional vials do not include any add-on products that regular-priced vials may include.

USE RESTRICTIONS

These cells are distributed for research purposes only.

This product is covered by a Limited Use License. By use of this product, the buyer agrees to the terms and conditions of all applicable Limited Use Label Licenses. For non-research use, such as screening, quality control or clinical development, contact outlicensing@invivogen.com.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: info@invivogen.com

