

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE  
1321N1**

---

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC0001-3

**Lot Number:** DC0001-3-061518

**Quantity:** 1 vial (4 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Amsbio Cellbanker 2 (11891)

**Host cell:** 1321N1

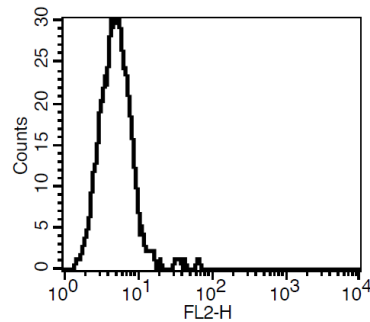
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS

**Stability:** Stable for 1-2 days after thawing

**Application:** Functional assays

**Figure 1**



**Figure 1.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thick line: parental cells.

**FOR RESEARCH USE ONLY.**

All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.