Primary Cells

Biologically Relevant Tools for Research

What Are They?

Non-transformed, non immortalized cells directly from the tissu



Why Use Them?

- Demonstrates tissue characteristics similar to in vivo conditions
- Emphasis from research community to use better cell culture tools
- Reduces animal testing costs if used for initial screens to refine experiments
- Limited lifespan maintains tissue-like characteristics versus cell lines that are passaged longer
- Valuable tool as qualified normal controls for primary diseased cell studies

Concerns with use of Cell lines

Authentication of cultured cell lines is critical for grant applications¹

are misidentified or cross-contaminated³

18-36% of cell lines

malignant prostatic cells and their normal epithelial counterparts are sought⁵

primary cultures of



500 leukemia-lymphoma cell lines, 15% of the cell lines were misidentified²

In a study of over

Some argue tumor cells obtained directly from patients are the best way to study cancer4

Nature News:

Primary Cells Limited lifespan retains cell identity Infinite Lifespan, loses cell specificity From the tissue From a vial with high mutations and clonal select 2





Mar M

- Your tools to get started on primary cell culture
 - Lonza's BulletKit™ MediumLonza's culturing protocols

Lonza's Primary Cells

- Lonza's scientific support team
- 1 Notice Number: NOT-0D-08-017 < grants.nih.gov/grants/guide/notice-files/NOT-0D-08-017.html>
- 2 Drexler HG: Guide to Leukemia-Lymphoma Cell Lines, 2nd Edition. Braunschweig, 2010, pages 883-892
 3 Hughes P et al. The costs of using unauthenticated, over-passaged cell lines: how much more data do we need? BioTechniques 43:575-586 (November 2007)
 4 Brendan Borrell. Nature News: How accurate are cancer cell lines?. Vol 463.18 February 2010.