

IncuCyte[®] MCF7 NuLight[™] Red

Catalog Number: 4524

Storage

Liquid Nitrogen

Note: Cells can be thawed and cultured immediately upon receipt or stored in liquid nitrogen for long-term storage. Storage at -80°C is not recommended. Cells should be used within 1 year of delivery.

Presentation

1mL, 1 x 10⁶ cells/mL in 90% FBS, 10% DMSO

Recommended Media and Components

- Eagle's Minimum Essential Medium (EMEM) (Cat# 30-2003 ATCC)
- 10% FBS (Cat# SH30071 Thermo Hyclone)
- 0.01mg/ml Human Recombinant Insulin (Cat# #12585014 Life Technologies)
- 1% Pen/Strep (Cat# 15140 Gibco/Life Technologies)
- 0.5µg/ml Puromycin (Cat# A11138-03 Gibco/Life Technologies)

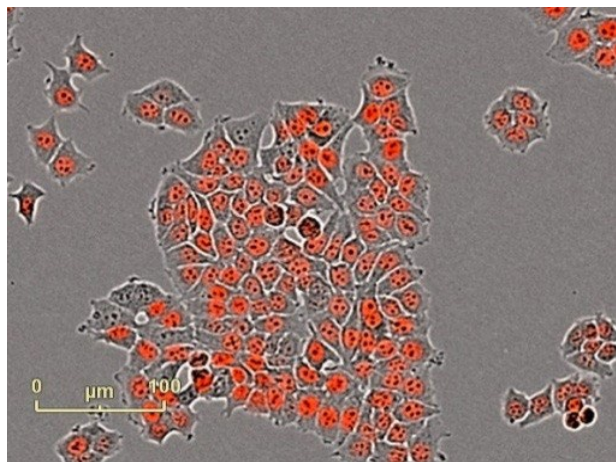
Background

Each vial contains a stable population of 1,000,000 MCF7 cells expressing the NuLight Red fluorescent protein restricted to the nucleus. Parental MCF7 cells were purchased from ATCC (Cat# HTB-22). MCF7 cells were transduced with the Essen NuLight Red Lentivirus (Cat# 4475; EF1α, puromycin) at an MOI of 3 (TU/cell) in the presence of 8µg/ml polybrene following the standard Essen protocol. This resulted in ≥70% transduction efficiency. 48 hours post infection, the complete population of cells was grown for 3-5 days in complete growth media containing 1µg/ml Puromycin to select for cells expressing NuLight Red.

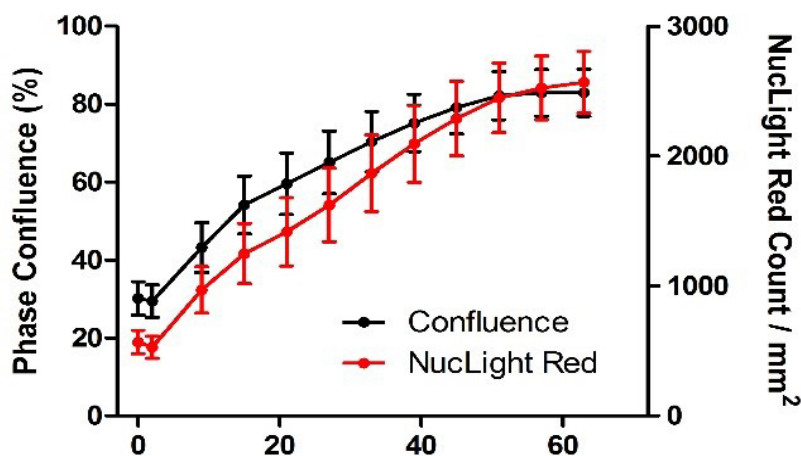
NuLight Red expressing cells are maintained in complete media containing 0.5µg/ml Puromycin. Following selection, a panel of validation assays designed to evaluate the effects of nuclear label expression on functional cell biology was completed. These assays include comparisons of cell morphology and growth/proliferation between stable populations and the parent populations from which they were derived (see below). In addition, all cells in our NuLight catalog have been certified mycoplasma free by ATCC and our stable populations have been unambiguously authenticated using ATCC's Short Tandem Repeat (STR) profiling.

Figure 1. MCF7 NuLight Red Cell Line

One vial of MCF7 NuLight Red cells was thawed into a 75cm² tissue culture treated flask and imaged in IncuCyte[®] ZOOM .



HD-phase contrast and red fluorescence blend.



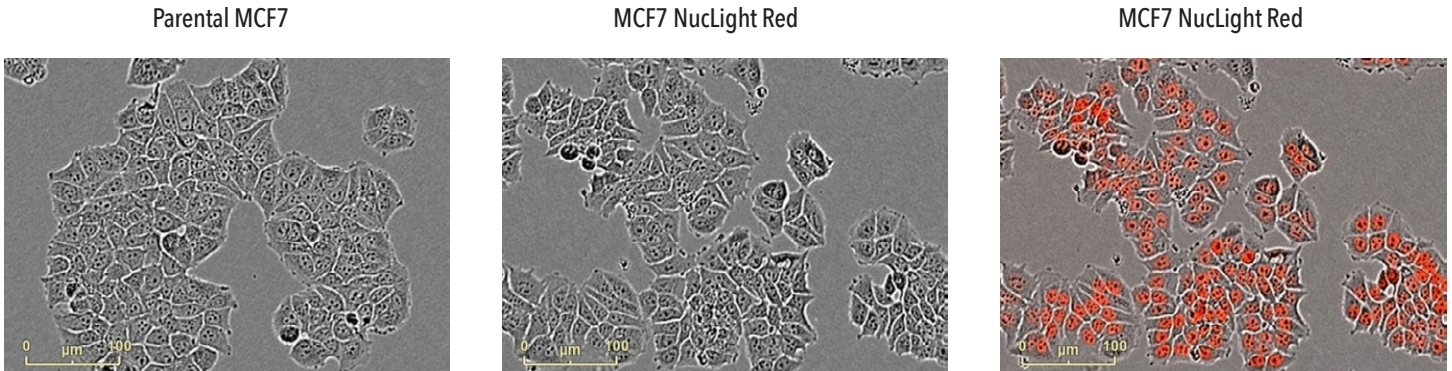
Phase confluence and nuclear counts over time.

MCF7 NuLight Red cells reach 80% confluence in a 75cm² flask after 2-3 days.

Validation Assays

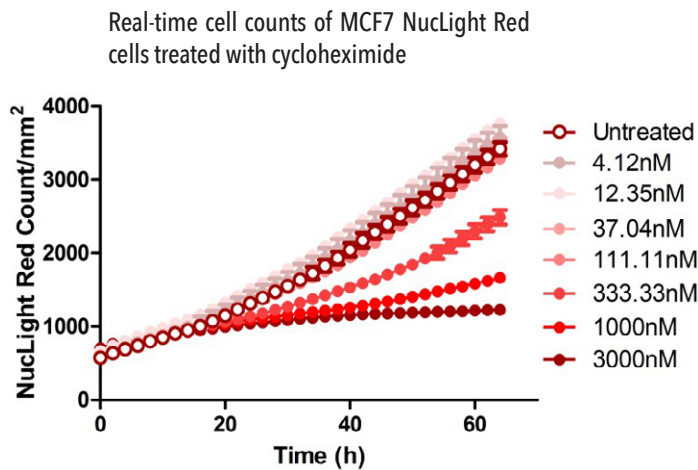
The following experiments were completed using an IncuCyte[®] ZOOM (10x)

1. Morphological comparison

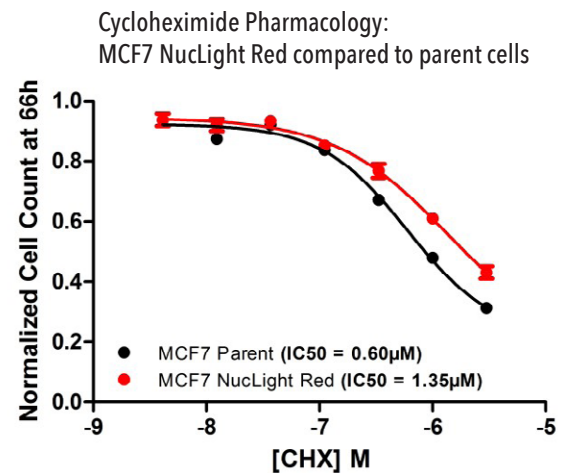


No significant alterations in cell morphology were observed between parental MCF7 cells and the MCF7 NuLight Red stable population.

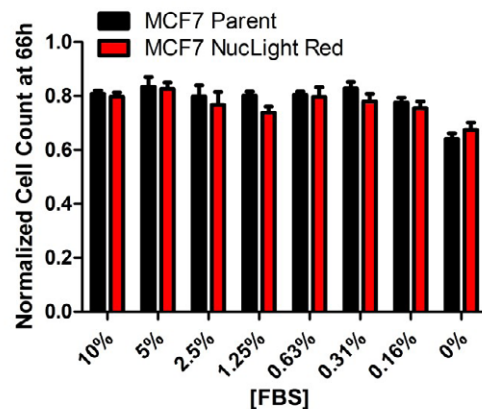
2. Proliferation – Real-time cell counts using the IncuCyte[®] Kinetic Proliferation Assay



Real-time cell counting revealed inhibition of MCF7 NuLight Red cell growth at cytostatic CHX concentrations ≥ 333 nM.



At the 48 hour endpoint, identically treated parental controls were stained with Vybrant DyeCycle Green and counted. Pharmacological analysis at the endpoint revealed similar CHX IC₅₀ concentrations for both parent and NuLight populations



Both Parent and NuLight populations were also grown in decreasing serum conditions. Endpoint values revealed similar growth characteristics in all concentrations.

Protocol:

Thawing and Culturing Cells

1. The recommended seeding density for MCF7 NuLight Red cells is 1,500 to 2,000 cells/cm². From this, calculate the number of flasks needed.
2. Prepare the flasks and culture media. To maintain expression of NuLight Red label, it is recommended that cells are maintained in complete media (EMEM + 0.01mg/ml Bovine Insulin + 10% FBS) containing 0.5µg/ml Puromycin. Puromycin can be removed for experiment/assay set-up.
3. Remove the vial containing MCF7 NuLight Red from liquid nitrogen.
4. In a 37°C water bath, quickly thaw the vial by gentle agitation. Be careful not to submerge entire vial to avoid contamination. This process should take no more than 2 minutes. Remove vial when only a tiny ice crystal remains.
5. Wipe vial with 70% ethanol. In a biosafety hood, aliquot cells into sufficient complete media to distribute among the flasks set up in Step 2.
6. Gently rock the culture flasks to evenly distribute the cells and place in 37°C incubator.

Related Products:

NuLight™/CytoLight™ Reagents

Cat.# 4475 CellPlayer NuLight Green (Lenti, EF-1 alpha, puro)

Cat.# 4481 CellPlayer CytoLight Green (Lenti, EF-1 alpha, puro)

Cat.# 4513 CellPlayer CytoLight Green (Lenti, CMV, no selection)

Cat.# 4476 CellPlayer NuLight Red (Lenti, EF-1 alpha, puro)

Cat.# 4482 CellPlayer CytoLight Red (Lenti, EF-1 alpha, puro)

NuLight Cell Lines:

Cat.# 4485 CellPlayer HT-1080 NuLight Red

Cat.# 4486 CellPlayer HT-1080 NuLight Green

Cat.# 4487 CellPlayer MDA-MB-231 NuLight Red

Cat.# 4488 CellPlayer MDA-MB-231 NuLight Green

Cat.# 4489 CellPlayer HeLa NuLight Red

Cat.# 4490 CellPlayer HeLa NuLight Green

Cat.# 4491 CellPlayer A549 NuLight Red

Cat.# 4492 CellPlayer A549 NuLight Green

Cat.# 4506 CellPlayer HUVEC NuLight Green

Cat.# 4511 CellPlayer Neuro-2a NuLight Green

Cat.# 4453 CellPlayer HUVEC CytoLight Green

Cat.# 4512 CellPlayer Neuro-2a NuLight Red

FOR RESEARCH USE ONLY. NOT FOR THERAPEUTIC OR DIAGNOSTIC USE.

Licences and warranty

Essen BioScience warrants that this product performs as described on the product label and in all literature associated with the sale of said product when used in accordance with the described protocol. This limited warranty is the sole warranty. No other warranties exist that extend beyond this warranty, either expressed or implied. Essen BioScience disclaims any implied warranty of merchantability or warranty of fitness for a particular purpose. Essen BioScience disclaims any responsibility for injury or damage and shall not be liable for any proximate, incidental or consequential damages in connection with this product.

If it is proven to the satisfaction of Essen BioScience that this product fails to meet performance specifications, Essen BioScience's sole obligation, at the option of Essen BioScience, is to replace the product or provide the purchaser with credit at or below the original purchase price. This limited warranty does not extend to anyone other than the purchaser. Notice of suboptimal performance must be made to Essen BioScience within 30 days of receipt of the product.

This Essen BioScience product contain proprietary nucleic acid(s) coding for proprietary fluorescent protein(s) being, including its derivatives or modifications, the subject of pending patent applications and/or patents owned by Evrogen JSC (hereinafter "Evrogen Fluorescent Proteins").

The purchase of Essen BioScience products incorporating these fluorescent proteins conveys to the buyer the non-transferable right to use Evrogen Fluorescent Proteins only for research conducted by the buyer. No rights are conveyed to modify or clone the gene encoding fluorescent protein contained in this product or to use Evrogen Fluorescent Proteins for commercial purposes. The right to use Evrogen Fluorescent Proteins specifically excludes the right to validate or screen compounds for commercial purposes. For information on commercial licensing, contact Evrogen Licensing Department, email: license@evrogen.com.