



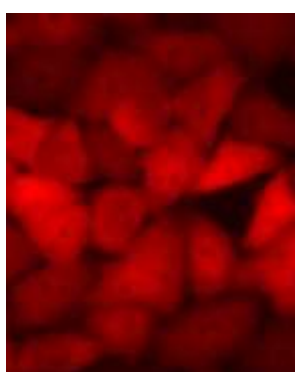
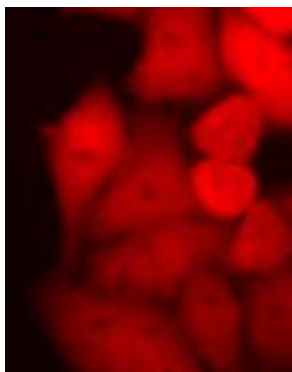
CellTrend GmbH

GESCHÄFTSFÜHRER: DR. HARALD HEIDECKE
IM BIOTECHNOLOGIEPARK • D-14943 LUCKENWALDE • E-MAIL: INFO@CELLTREND.DE
TEL: +49 (0)3371 / 681 290 • FAX: +49 (0)3371 / 681 312

Cell Line Specification Sheet

HeLa-TurboFP602

Human cervix carcinoma cells expressing deep-red fluorescent TurboFP602

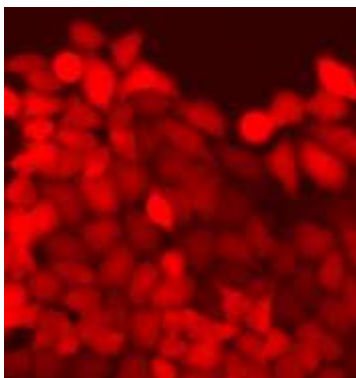


SPECIFICATIONS:

Cell Line:	HeLa cells expressing TurboFP602
Cells/vial:	approximately 5×10^5
Subculturing:	1:3 to 1:5
Medium Renewal:	2 to 3 times per week
Growth Medium:	DMEM (1 g/L D-Glucose, 3,7 g/L NaHCO ₃) + 10 % fetal calf serum (FCS), 2 mM L-Alanyl-L-Glutamine, 1 % Non Essential Amino Acids (NEA)
Freeze Medium:	complete growth medium + 10% FCS, 10% DMSO
mycoplasma-free	
Shipping:	frozen
Storage Recommendation:	Liquid nitrogen
Biosafety Level:	1

INFORMATION:

Organism:	<i>homo sapiens</i> (human)
Source:	Organ: cervix, disease: adenocarcinoma
Gender:	female, 31 years
Ethnicity:	African American
Growth properties:	adherent
Morphology:	epithelial
Conditions:	37 °C, 5 % CO ₂
Plasmid:	pTurboFP602-C (Evrogen, Moscow, Russia)



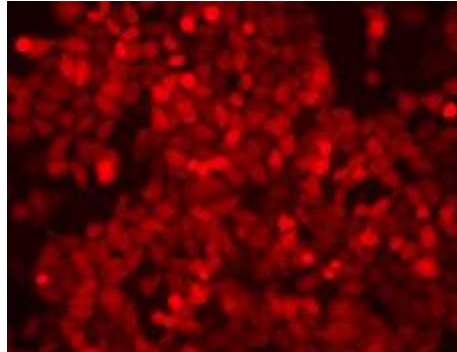
HeLa cells expressing TurboFP602

Human cervix carcinoma cells were stably transfected with pTurboFP602 resulting in a bright red fluorescence with an emission maximum at 602 nm in all cells. The HeLa cell line was established as the first cell line from a cervix carcinoma of a 31-year-old woman in 1951. These cells proliferate rapidly and display an active telomerase enzyme avoiding normal aging and cell death through shortening of the telomeres. HeLa cells are very durable, easy to handle and widely used in cell culture-based research.

FOR NON-HUMAN INVESTIGATIONAL RESEARCH ONLY.

TO BE HANDLED UNDER BIOSAFETY LEVEL 1 CONTAINMENT.

Information derived from ATCC and ECACC.



This product contains a proprietary nucleic acid from coding for a proprietary fluorescent protein(s) intended to be used for research purpose only, where research purpose means uses or activities (or the results therefrom) that do not directly generate revenue. Any use of the proprietary nucleic acid or protein other than for research use is strictly prohibited. USE IN ANY OTHER APPLICATION REQUIRES A LICENSE FROM EVROGEN. To obtain such a license, please contact Evrogen at license@evrogen.com.

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com • info@eaglebio.com



EAGLE
BIOSCIENCES