



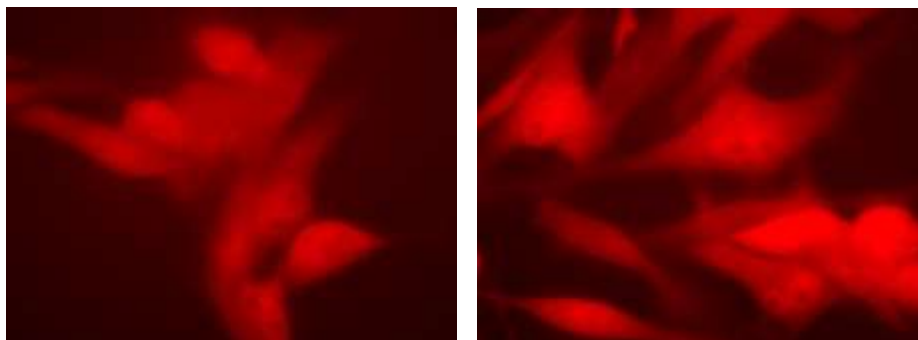
CellTrend GmbH

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

M3-JRed

Mouse melanoma cells expressing JRed

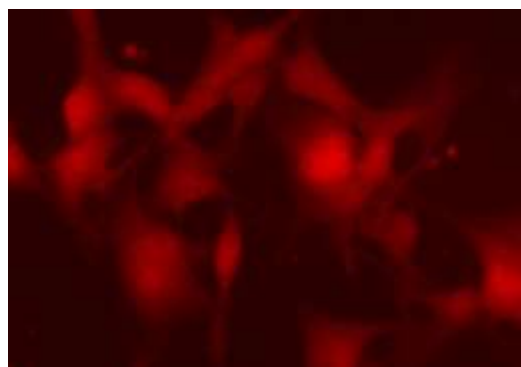


SPECIFICATIONS:

Cell Line:	M3 cells (clone of Cloudman S91-M3 melanoma) expressing the red fluorescent protein JRed
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>Mus musculus</i> (mouse)
Source:	Organ: skin; Disease: melanoma
Gender:	male
Tumorigenic:	yes
Growth properties:	adherent
Morphology:	epithelial
Conditions:	37 °C, 5 % CO ₂
Plasmid:	pJRed (Evrogen, Moscow, Russia)



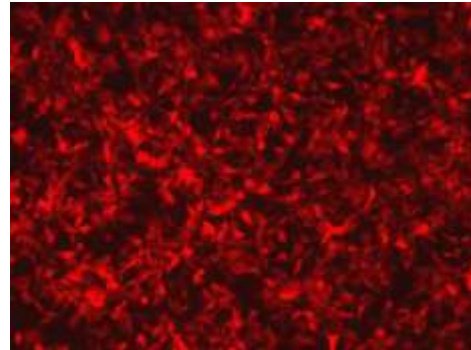
M3 cells expressing JRed

A bright red fluorescent cell line has been established by stable transfection with a plasmid coding for JRed. The M3 mouse melanoma (clone of Cloudman S91-M3) cell line constitutes an ideal model for in vivo experiments. Melanoma-like tumors in different colors can be induced in syngeneic mice by subcutaneous injection of M3-JRed and M3-PhiYFP cells. After several days tumor formation becomes evident.

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