

Customer recommended DharmaFECT™ transfection reagent formulation for various cell lines

The information provided in the table below represents anecdotal customer feedback from researchers reporting successful siRNA transfection with one or more DharmaFECT™ transfection reagent formulations. Please note that our scientists have not validated this

Human		
Cell line	Cell type	Successful DharmaFECT formulation
143B	Osteosarcoma	1
786-0	Renal epithelial adenocarcinoma	4
A-431	Skin Epidermal Carcinoma	1
AU565	Breast Adenocarcinoma	1
BT-474	Breast Carcinoma	1
Caco-2	Colorectal Adenocarcinoma	1
DLD-1	Colorectal Adenocarcinoma	4
GM5756T	Fibroblasts	4
H1155	Lung Carcinoma	1
H2122	Lung Adenocarcinoma	1, 2, 3, 4
H2126	Lung Adenocarcinoma	3
hASMC	Aortic Smooth Muscle	1
HBEC30	Lung Carcinoma	3
HBEC34	Lung Carcinoma	2
HBEC3-KT	Lung Carcinoma	1
HCC1569	Breast Carcinoma	1
HCC193	Nonsmall cell lung carcinoma	1, 2, 3, 4
HCC1954	Breast Carcinoma	1
HCC366	Lung Adenocarcinoma	1

information and offers this as a courtesy. We encourage customers to optimize important parameters such as cell number, DharmaFECT volume, and control siRNA. See the [DharmaFECT General Transfection Protocol](#) for optimization guidelines.

Human		
Cell line	Cell type	Successful DharmaFECT formulation
HCC4017	Lung Carcinoma	2
HCT116	Colorectal Carcinoma	1, 2, 3, 4
Hep3B	Hepatocellular Carcinoma	4
HEY	Ovarian Carcinoma	1
HL60	Promyelocytic Leukemia	1
Hs578T	Breast Adenocarcinoma	1
Huh7	Hepatoma	1, 4
IGROV1	Ovarian Carcinoma	1
IHGK	Immortal Keratinocytes	1
IMR-32	Neuroblastoma	2
Jurkat	Leukemia T-cell lymphoblast	4
LS180	Colon Carcinoma	1
MDA-MB 361	Breast Adenocarcinoma	1
MDA-MB-231	Breast Adenocarcinoma	1
MIA PaCa-2	Pancreatic Carcinoma	2
MS53	Lung Carcinoma	4
OVCAR-3	Ovarian Carcinoma	1
OVCAR-5	Ovarian Carcinoma	1
OVCAR-8	Ovarian Carcinoma	1

Human		
Cell line	Cell type	Successful DharmaFECT formulation
PANC-1	Pancreatic Carcinoma	4
PANC-10	Pancreatic Carcinoma	4
Primary	Primary Chondrocytes	1, 2, 3, 4
Primary	Primary Fibroblasts	1, 3
Primary	Primary Microglial	3
Primary	Primary Placental Trophoblasts	2
SCLC	Small Cell Lung Carcinoma	2
SKBR3	Breast Epithelial Adenocarcinoma	1
SK-OV-3	Ovarian Adenocarcinoma	1, 2, 3, 4
SKW6.4	B lymphocyte	4
SNB19	Glioblastoma	4
SW 1353	Chondrosarcoma	1
SW620	Colorectal Adenocarcinoma	1, 2, 3
T98G	Gioblastoma	2
U-2 OS	Osteosarcoma	1
U251	Glioblastoma	4
UACC816	Breast carcinoma	1
WI-38	Lung fibroblast	1
ZR-75-30	Breast carcinoma	1

Mouse		
Cell line	Cell type	Successful DharmaFECT formulation
3T3-Swiss	Fibroblast	3
COMMA-D	Mammary Epithelial	1
G1E	Gata-1 erythroid-committed cell line	2
MEF	Embryonic Fibroblast	4
Neuro2A	Neuroblastoma	2
NIH/3T3	Embryonic Fibroblast	3
NMuMG cells	Mammary Epithelial	4
Primary	Primary Glial	3
Primary	Primary Muscle	4
Primary	Primary Hepatocytes	1
Primary	Primary Neuronal	3
RGC	Retinal Ganglion	4
S91	Cloudman S91 Melanoma	1

Rat		
Cell line	Cell type	Successful DharmaFECT formulation
Cardiac myocytes	Neonatal - Primary	1
FaO	Liver Hepatoma	2
H9c2	Cardiomyoblast	1,2
L2	Lung Epithelial	1
L6	Skeletal Muscle Myoblast	3
McA-RH7777	Liver Hepatoma	4
PC12	Adrenal Gland Pheochromocytoma	1, 2, 3
Primary	Primary Cardiac Myocytes	1
Primary	Primary Hippocampal	3
Primary	Primary Hepatocytes	1
Primary	Primary Neuronal	3
RC4B	Lactogonadotrope	1, 2, 3, 4

Other		
Cell line	Cell type	Successful DharmaFECT formulation
Primary	Canine Primary chondrocyte	2,4
Primary	Bovine Microvascular Retinal Endothelial	1
LR73	Chinese Hamster Ovary Derivative	2

If you have any questions, contact

t +44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA)

f +44 (0)1223 655 581

w horizondiscovery.com/contact-us **or** dharmacon.horizondiscovery.com/service-and-support

Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

All trademarks are the property of Horizon Discovery Company unless otherwise specified. ©2018 Horizon Discovery Group Company—All rights reserved. First published January 2015. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.

horizonTM
INSPIRED **CELL** SOLUTIONS

V20818