

PART NO.	PRODUCT DESCRIPTION	UNIT SIZE	UNIT PRICE
<b>HUMAN CELLS AND MEDIA</b>			
NHEK-CRY-NEO	Normal Human Epidermal Keratinocytes - Neonatal, Cryopreserved	>500K Cells	\$586.00
NHEK-CRY-AD	Normal Human Epidermal Keratinocytes - Adult, Cryopreserved	>500K Cells	\$655.00
NHEK-GM	Normal Human Epidermal Keratinocytes - Growth Medium	500 ml	\$148.00
NHEK-BM	Normal Human Epidermal Keratinocytes - Basal Medium	500 ml	\$99.00
NHEK-GS	Normal Human Epidermal Keratinocytes - Growth Supplement	5 ml	\$77.00
NHM-CRY-NEO	Normal Human Melanocytes - Neonatal, Cryopreserved	>500K Cells	\$792.00
NHM-CRY-AD	Normal Human Melanocytes - Adult, Cryopreserved	>500K Cells	\$655.00
NHM-BM	Normal Human Melanocytes - Basal Medium	500 ml	\$99.00
NHM-GM	Normal Human Melanocytes - Growth Medium	500 ml	\$153.00
NHM-GS	Normal Human Melanocytes - Growth Supplement	10 ml	\$105.00
NHDF-CRY-NEO	Normal Human Dermal Fibroblasts - Neonatal, Cryopreserved	>500K Cells	\$655.00
NHDF-CRY-AD	Normal Human Dermal Fibroblasts - Adult, Cryopreserved	>500K Cells	\$449.00
NHDF-BM	Normal Human Dermal Fibroblasts - Basal Medium	500 ml	\$119.00
NHDF-GM	Normal Human Dermal Fibroblasts - Growth Medium	500 ml	\$153.00
NHDF-GS	Normal Human Dermal Fibroblasts - Growth Supplement	5 ml	\$77.00
<b>HUMAN AIRWAY CELLS</b>			
NHBE-CRY	Normal Human Bronchial Epithelial Cells, Cryopreserved	>500K Cells	\$959.00
NHBE-CRY-DS	Human Bronchial Epithelial Cells - Diseased (Asthma, COPD, Smoker)	>500K Cells	\$1,136.00
NHBE-BM	Normal Human Bronchial Epithelial Cells - Basal Medium	500 ml	\$119.00
NHBE-GM	Normal Human Bronchial Epithelial Cells - Growth Medium	500 ml	\$182.00
NHBE-GS	Normal Human Bronchial Epithelial Cells - Growth Supplement	10 ml	\$112.00
NHBE-HCS	Normal Human Bronchial Epithelial Cells - Aliquot for supplementation of NHBE-BM	250 µl	\$20.00
NHBF-CRY	Normal Bronchial Fibroblast Cells, Cryopreserved	>500K Cells	\$758.00
NHBF-CRY-DS	Human Bronchial Fibroblast Cells - Diseased (Asthma, COPD, Smoker)	>500K Cells	\$861.00
<b>HUMAN OCULAR CELLS</b>			
NHCE-BM	Normal Human Corneal Epithelial Cells - Basal Medium	500 ml	\$104.00
NHCE-CM	Normal Human Corneal Epithelial Cells - Coating Mixture	10 ml	\$70.00
NHCE-CRY	Normal Human Corneal Epithelial Cells, Cryopreserved	>500K Cells	\$861.00
NHCE-GM	Normal Human Corneal Epithelial Cells - Growth Medium	500 ml	\$137.00
NHCE-GS	Normal Human Corneal Epithelial Cells - Growth Supplement	5 ml	\$77.00
NHCK-CRY	Normal Human Corneal Keratocytes (Corneal Fibroblasts) - Cryopreserved	>500K Cells	\$861.00

www.mattek.com | support@mattek.com

**USA Headquarters**  
200 Homer Ave  
Ashland, MA 01721

T: +1-508-881-6771  
F: +1-508-879-1532

PART NO.	PRODUCT DESCRIPTION	UNIT SIZE	UNIT PRICE
<b>HUMAN IMMUNE CELLS</b>			
DC-100-CRY	Human Dendritic Cells - Cryopreserved	>2.5M Cells	\$930.00
DC-100-RT	Human Dendritic Cells - Ready To Use	>2.5M Cells	\$993.00
DC-BAS-50	Human Dendritic Cells - Basal Medium, 50 ml	50 ml	\$99.00
DC-BAS-250	Human Dendritic Cells - Basal Medium, 250 ml	250 ml	\$236.00
DC-MM-50	Human Dendritic Cells - Maintenance Medium, 50 ml	50 ml	\$148.00
DC-MM-250	Human Dendritic Cells - Maintenance Medium, 250 ml	250 ml	\$663.00
DCP-100-CRY	Human Plasmacytoid Dendritic Cells - Cryopreserved	>2.5M Cells	\$1,081.00
DCP-100-RT	Human Plasmacytoid Dendritic Cells - Ready To Use	>2.5M Cells	\$1,158.00
DCP-MM-50	Human Plasmacytoid Dendritic Cells - Maintenance Medium, 50 ml	50 ml	\$208.00
DCP-MM-250	Human Plasmacytoid Dendritic Cells - Maintenance Medium, 250 ml	250 ml	\$765.00
TC-100-CRY	Human T-Cells - Cryopreserved	>2.5M Cells	\$993.00
TC-100-RT	Human T-Cells - Ready To Use	>2.5M Cells	\$1,041.00
TC-BAS-50	Human T-Cells - Basal Medium, 50 ml	50 ml	\$99.00
TC-BAS-250	Human T-Cells - Basal Medium, 250 ml	250 ml	\$236.00
TC-MM-250	Human T-Cells - Maintenance Medium, 250 ml	250 ml	\$765.00

[www.mattek.com](http://www.mattek.com) | [support@mattek.com](mailto:support@mattek.com)

**USA Headquarters**  
 200 Homer Ave  
 Ashland, MA 01721

T: +1-508-881-6771  
 F: +1-508-879-1532