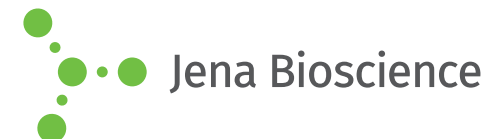




JBScreen Plus Volatiles

Cat.-No.: CS-505

SCREEN FORMULATION



No.	Additive	Type	CAS Number
V/A1	40.5 % v/v Methanol	Small Alcohol	67-56-1
V/A2	29.2 % v/v Ethanol	Small Alcohol	64-17-5
V/A3	38.2 % v/v 2-Propanol	Small Alcohol	67-63-0
V/A4	37.4 % v/v 1-Propanol	Small Alcohol	71-23-8
V/A5	8.2 % v/v 1-Butanol	Small Alcohol	71-36-3
V/A6	47.8 % v/v 2-Methyl-2-propanol	Small Alcohol	75-65-0
V/B1	27.9 % v/v Ethylene glycol	Small Alcohol	107-21-1
V/B2	36.1 % v/v 1,3-Propanediol	Small Alcohol	504-63-2
V/B3	35.4 % v/v 1,4-Butanediol	Small Alcohol	110-63-4
V/B4	35.8 % v/v 1,2-Butanediol	Small Alcohol	584-03-2
V/B5	29.1 % v/v 2,2,2-Trifluoroethanol	Small Alcohol	75-89-8
V/B6	21.1 % v/v 1,1,1,3,3,3-Hexafluoro-2-propanol	Small Alcohol	920-66-1
V/C1	30.8 % v/v 4-Butyrolactone	Non ionic Chaotrope	96-48-0
V/C2	52.2 % v/v Acetonitrile	Non ionic Chaotrope	75-05-8
V/C3	10.7 % v/v Propionitrile	Non ionic Chaotrope	107-12-0
V/C4	4.8 % v/v Methyl tert-butyl ether	Non ionic Chaotrope	1634-04-4
V/C5	34.1 % v/v 1,4-Dioxane	Non ionic Chaotrope	123-91-1
V/C6	32.4 % v/v Tetrahydrofuran	Non ionic Chaotrope	109-99-9
V/D1	36.7 % v/v Acetone	Organic Volatile	67-64-1
V/D2	22.4 % v/v Butanone	Organic Volatile	78-93-3
V/D3	3.7 % v/v 3-Pentanone	Organic Volatile	96-22-0
V/D4	4.00 % v/v Methyl acetate	Organic Volatile	79-20-9
V/D5	2.9 % v/v Ethyl acetate	Organic Volatile	141-78-6
V/D6	0.3 % v/v Dichloromethane	Organic Volatile	75-09-2

