

CloneMatrix

Methylcellulose concentrate for semi-solid media

KEY FEATURES

- Add directly to already optimized media
- Animal-free
- Conveniently provided as 40 ml of 2.5X concentrated product in a 100 ml bottle

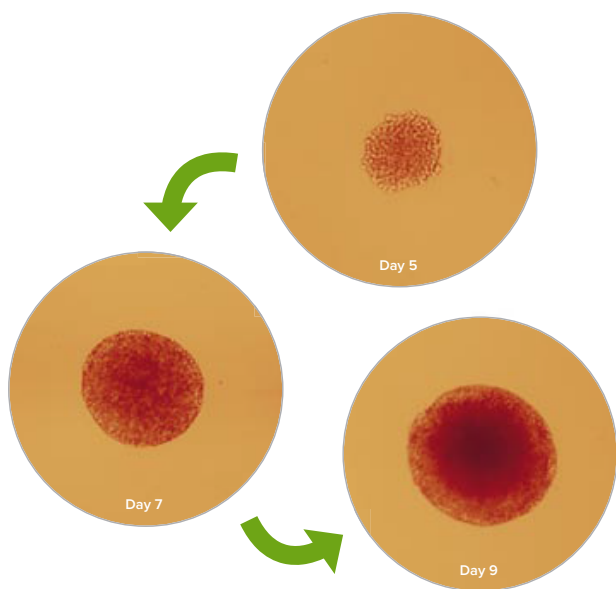
CloneMatrix™ is an easy-to-use product designed to convert existing mammalian cell biology media to semi-solid form. Semi-solid media serves to maintain the positional integrity of growing cell colonies thus allowing the automated identification and picking of clonal cell lines.

The use of semi-solid media combined with the unique fluorescent screening and picking technology of the ClonePix System will revolutionize clonal cell line development in biopharmaceuticals.

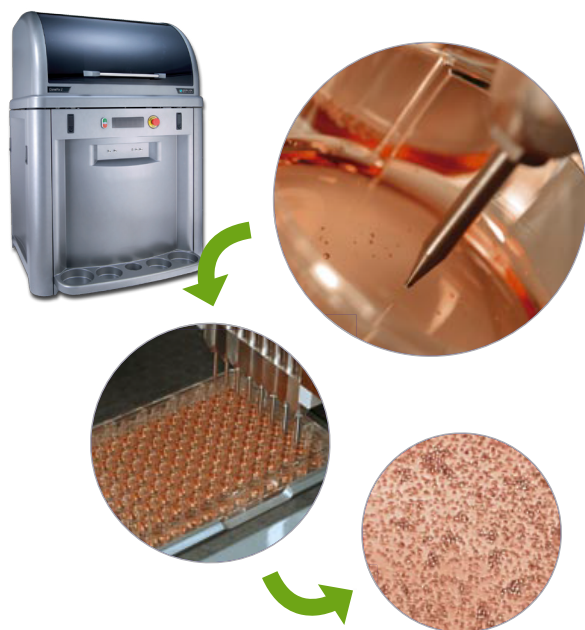
- CloneMatrix is a unique methylcellulose concentrate which offers the user flexibility to add their own media formulation to produce an optimized semisolid media reagent.
- CloneMatrix is conveniently supplied as 40 mL in a 100 mL bottle, reducing the risk of contamination as there is no requirement to aliquot the concentrate. Supplied as a 2.5x concentrate, simply add optimized media (2x concentration), selection and supplements as required.
- CloneMatrix is supplied as an animal-free product, so is suitable for use with serum-free media.

Ordering information

Description	Volume	Cat no.
CloneMatrix Methylcellulose concentrate for semi-solid media	40 mL (for 100 mL media)	K8510
CloneMatrix Methylcellulose concentrate for semi-solid media	6 x 40 mL (each for 100 mL media)	K8500
CloneMatrix optimized for CHO cells	40 mL (for 100 mL media)	K8530
CloneMatrix optimized for CHO cells	6 x 40 mL (each for 100 mL media)	K8520



The images show NS1 hybridoma colonies growing in CloneMatrix containing media. The images were taken 5, 7 and 9 days post seeding respectively. Images are at 40x magnification.



CloneMatrix media is optimized to work with ClonePix Systems.

Contact Us

Phone: [+1-800-635-5577](tel:+18006355577)
 Web: www.moleculardevices.com
 Email: info@moldev.com
 Check our website for a current listing of worldwide distributors.