



Optimize Screening for Biotherapeutics

DiSH™ Kit (ENZ-71001)

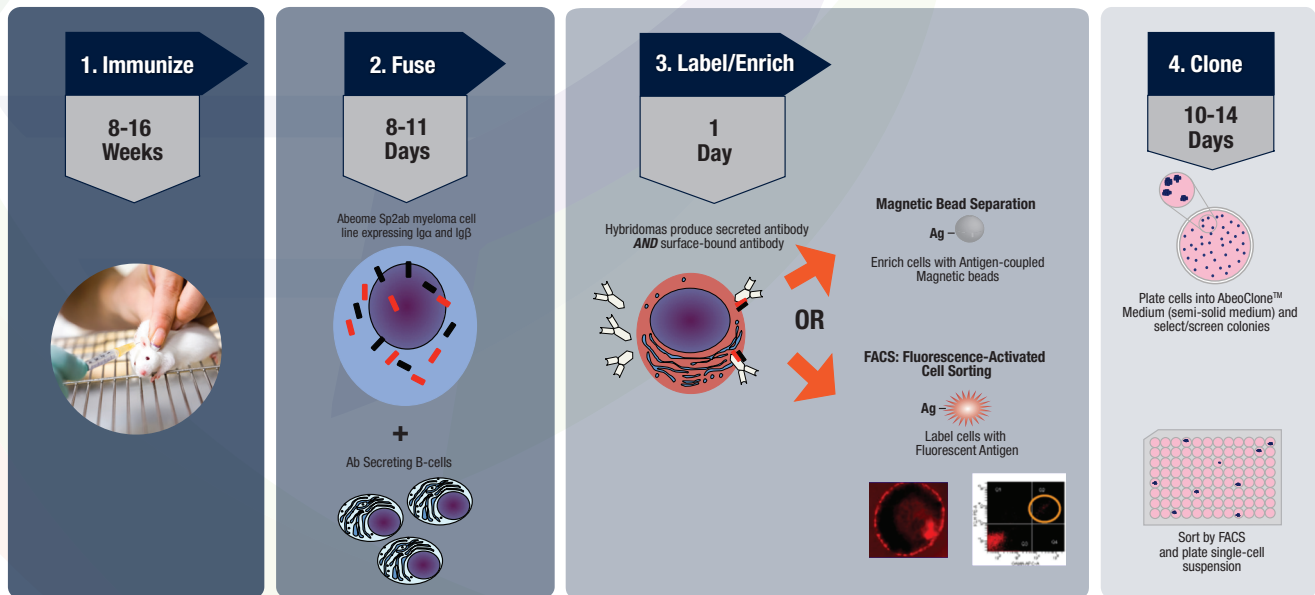
Reduce labor time and costs using a novel technology for rapid hybridoma production

The Direct Selection of Hybridomas (DiSH™) kit is a novel system in which mAb-secreting hybridoma cells co-express significant amounts of the membrane form of the secreted immunoglobulin (Ig) on their surfaces. Subsequently, hybridomas with high-affinity for the target antigen are labeled. Once labeled, these hybridomas can be isolated from among a pool of hundreds of thousands of cells in a matter of hours using Fluorescence Activated Cell Sorting (FACS) or magnetic separation. The total elapsed time from harvest to the first screening is typically 16 days, resulting in improved efficiency for detection of highly specific mAb.

Key Features

- Unique technology for more rapid hybridoma production
- Decreased time and labor costs
- Improved selection of specific, stable hybridomas
- Minimal manipulation of cloned cell lines when compared to approaches requiring limiting dilution subcloning (LDS)

PRODUCT NAME	PRODUCT #
DiSH™ Kit	ENZ-71001
AbeoClone™ (Base Semi-solid Medium)	ENZ-70009
AbeoClone™ (HAT Semi-solid Medium)	ENZ-70004
AbeoClone™ (CHO TCSC Semi-solid Medium)	ENZ-70010
Sp2ab Myeloma Fusion Partner	ENZ-70008



Abeome products and product applications are covered by US Patent No. 7,148,040.

Accelerate Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com