

## CLONEPIX™ 2 APPLICATIONS TIPS & RECOMMENDATIONS FOR SUCCESSFUL PICKING

### Prior to Picking:

#### Plates & Media:

- Ensure 2mL of media per well in 6-well plate, with media touching edges of wells. Do not disturb plated cells for at least 5 days post plating
  - Disposable pipettes may be useful in more accurate dispensing.
- Keep well plates level to ensure even distribution of media during incubation
  - Different media heights can affect imaging readings
- Keep plates steady during transport
- Add sterile water to reservoirs around wells to retain humidity and prevent the evaporation of the semisolid media.

#### Wash Bottles:

- Ensure all wash bottles are full and caps are secure
- To reduce contamination, use autoclaved/sterilized water for picking.
  - Millex-LG 0.20µm filters (included) help minimize sterility issues.
- When changing wash bottles, run the “Sanitize Pins” function under Utilities on the Fusion Software Main Page for a minimum of 30 cycles for the head to purge

Sanitize Options

Number of cycles for the head to purge:

Number of cycles the pins will be scrubbed in the Wash Bath:

Number of seconds the Dryer will be on for:

#### Picking Head:

- F2 (0.7mm) Pins are recommended for Adherent Cell Picking
- F1 (0.4mm) Pins are recommended for Suspension Cell Picking
- When changing heads or pins, run the “Picking Head Alignment” function under Utilities on the Fusion Software Main Page

#### Picking Settings:

Pin Options:

Adherent

Suspension

Normal Aspirate/Dispense Volumes

Aspirate Volume – 5 to 8 µL

Dispense Volume – 7 to 10 µL  
(Equal to Aspirate volume + 2 µL)

Semi-Solid Media Delay (ms):

Pick Height Adjustment:  Above Well Bottom

Aspirate Volume (µl):

Dispense Volume (µl):

Sanitize Options

**Recommended Wash Settings**

Number of cycles for the head to purge:

Number of cycles the pins will be scrubbed in the Wash Bath:

Number of seconds the Dryer will be on for:

#### Group Settings:

Feature	F1 Pins	F2 Pins
Too Big	0.7	0.7
Too Small	0.03	0.03
Irregular 1	0.6	0.6
Irregular 2	0.6	0.6
Proximity*	0.7-1.0	1.0-1.2

\*For better monoclonality assurance, a minimum of 1.0mm is recommended.

#### And Always Remember...

