

Update: SOLiD™ ToP Paired End 50/35 bp Sequencing Kits

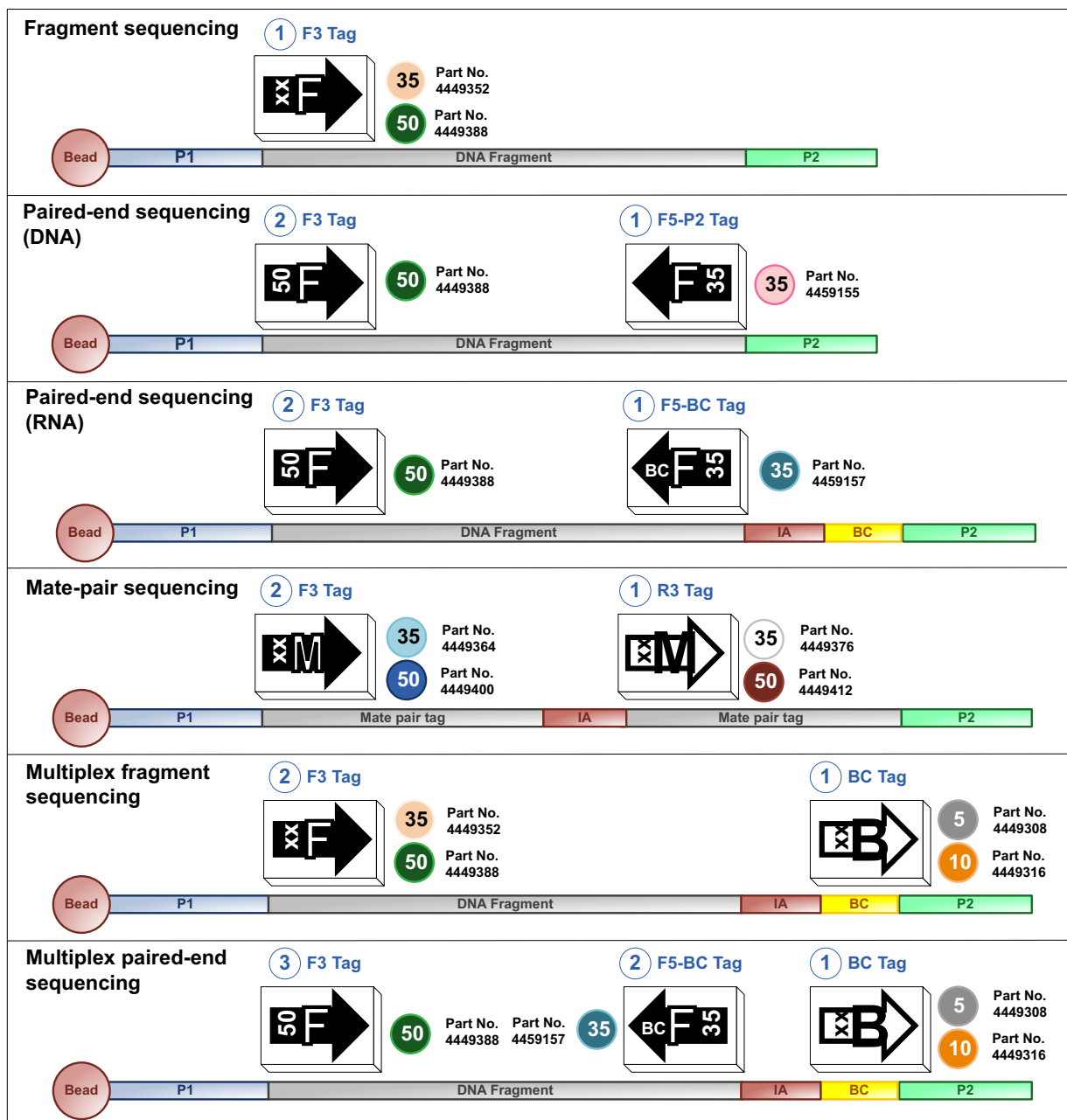
Note: For safety and biohazard guidelines, refer to the “Safety” section in the *Applied Biosystems SOLiD™ 4 System Instrument Operation Guide* (PN 4448379). For every chemical, read the SDS and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

The updated kit loading order and reagent strip layout should be used with the new 50/35-bp paired-end sequencing kits.

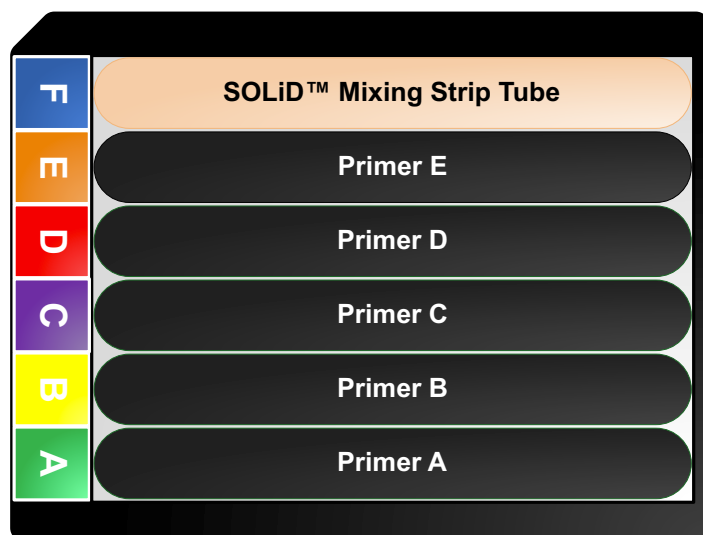
Kit loading order

Reagent strip box symbols correspond with sequencing of fragment, mate-paired, and barcoded fragment libraries. For more information, refer to the *Applied Biosystems SOLiD™ 4 System Product Selection Guide* (PN 4452360).

Note: Verify that there is sufficient volume of Reset Buffer for each tag sequenced.



Paired-end sequencing (F5-P2/F5-BC tags) reagent strip layout



SOLiD™ ToP Paired End Seq. Kit – Frag. Lib., F5-P2 Tag, MM35

SOLiD™ ToP Paired End Seq. Kit – BC Frag. Lib., F5-BC Tag, MM35

For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.

NOTICE TO PURCHASER: PLEASE REFER TO THE APPLIED BIOSYSTEMS SOLiD™ 4 SYSTEM INSTRUMENT OPERATION GUIDE FOR LIMITED LABEL LICENSE OR DISCLAIMER INFORMATION.

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

© 2010 Life Technologies Corporation. All rights reserved.

Part Number 4458612 Rev. A 08/2010

