

## More answers from a single tub

BD Multitest™ 6-color TBNK reagent is indicated for use in the immunological assessment of normal individuals, and patients having, or suspected of having, immune deficiency. BD Multitest™ 6-color TBNK with optional BD Trucount™ tubes is intended for use with BD FACSLyric™, BD FACSCanto™ II, and BD FACSCanto™ flow cytometers to determine the percentages and absolute counts of mature human lymphocyte subsets in peripheral whole blood for immunophenotyping.

