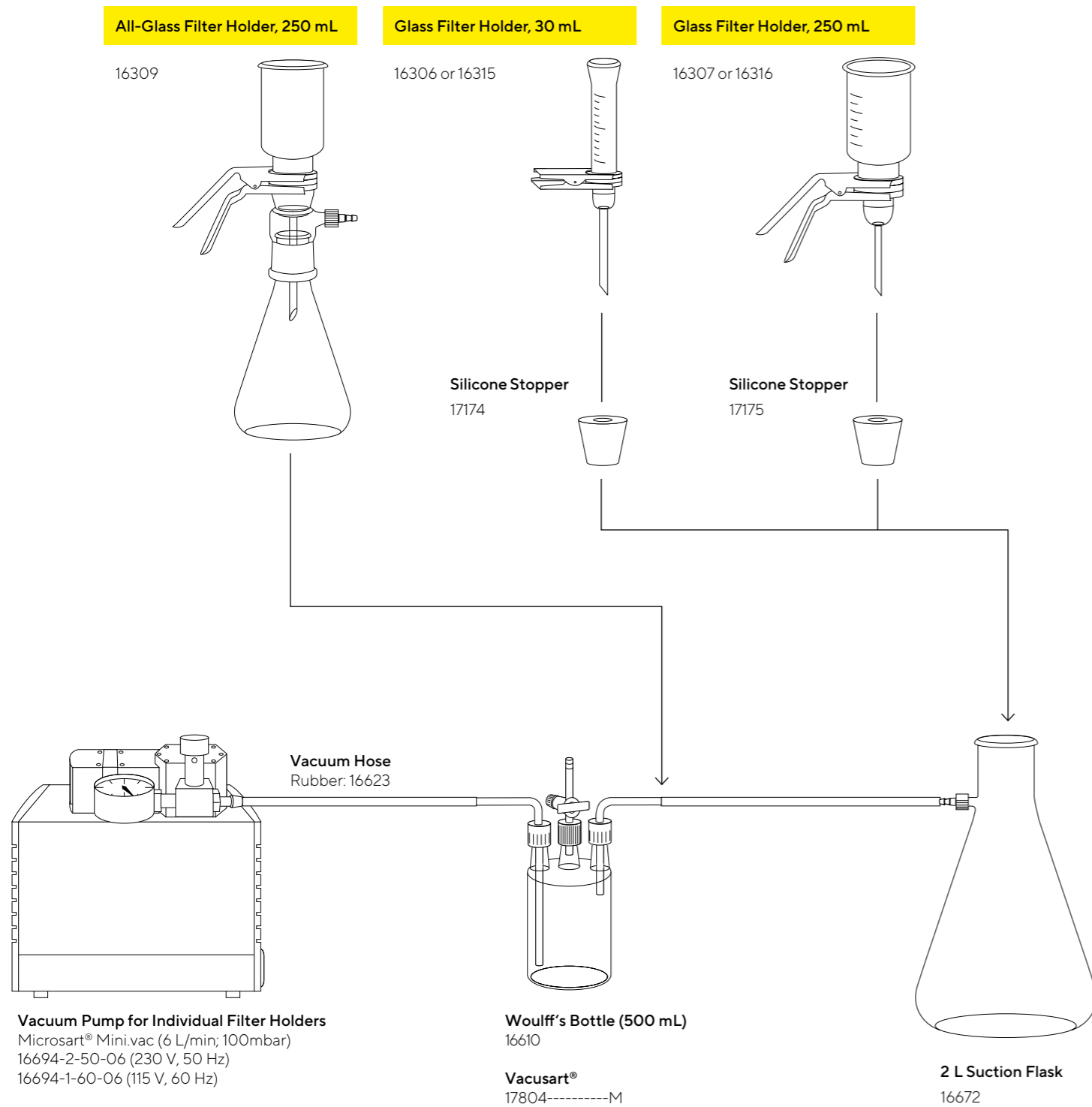


Selection of Filtration Accessories for Parallel Vacuum Filtration

Individual Filter Holders



3-Branch Manifold

