

➤ Field of Application

Clinical data indicate that only a part of patients respond to treatment with immune-modulating drugs. It has therefore been recognized as important to consider the immune response already during preclinical drug development. Flow cytometry is widely used for analysis of immune cell populations in tumors.

➤ All-In-One Flow Cytometry Panel (17 markers)

A comprehensive analysis of numerous immune cell populations in one tumor is a major challenge. We use one staining panel that allows for the analysis of all major immune cell populations and comprises live/dead dye and antibodies against CD3, CD4, CD8a, CD45, CD25, CD11b, Ly6C, Ly6G, F4/80, CD11c, MHC class II, CD206, CD335, CD49b, B220 and FoxP3.

