



NOTE: This poster conveys a general overview and should be considered neither comprehensive nor definitive. The details of this information are understood to be subject to interpretation. © R&D Systems 2016

Learn more | rndsystems.com/microgliaposter