

PRE-PRIMING
(for non-canonical inflammasomes)

PRIMING (signal 1)

ACTIVATION (signal 2)

OUTCOMES

OVERVIEW OF INFLAMMASOME SENSORS AND ACTIVATORS

