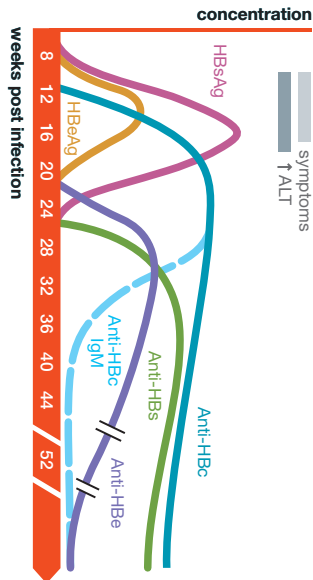
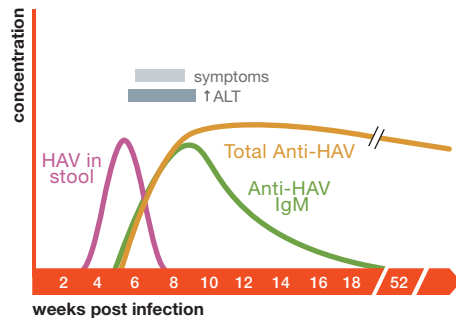


Hepatitis Serology

Hepatitis B



Hepatitis A



Hepatitis C

