濃度がずれる場合のチェックポイント







ピークがスプリットしていたり、位置がずれている場合には



a)グラフ上で右クリック>Manual Integration選択

b)右クリック>Add Peak, Remove Peakにてピーク追加・削除

c)青線で各ピークごとのベースライ ンが表示されるようになるので、 ベースライン末端の終点・始点を ドラッグで動かす



ソフトウェアがUpperMarkerピークを正しく捉えていない場合





正しいピークを選択し、右ク リック>Manually Set Upper Markerを選択して下さい。

